



Eddy Covariance Publications List

Citations related to the use of LI-COR gas analyzers and software in eddy covariance experiments

This document contains a sampling of recent publications that reference LI-COR instrumentation and software. This list is provided for informational purposes only, and LI-COR neither endorses, nor makes any express or implied warranties with respect to any data included in these publications.

1 January 2014 to 1 October 2015

Volume 1

LI-7700 Open Path CH₄ Analyzer — below
LI-7200 Enclosed CO₂/H₂O Analyzer — page 2
LI-7500/A Open Path CH₄ Analyzer — page 5
EddyPro® Eddy Covariance Processing Software — page 20

LI-7700 Open Path CH₄ Analyzer

- Alberto, M. C. R., et al. (2014). **Measuring methane flux from irrigated rice fields by eddy covariance method using open-path gas analyzer.** *Field Crops Research* 160: 12–21.
- Begashaw, I., et al. (2014). **A new tool for automated data collection and complete on-site flux data processing for eddy covariance measurements.** AGU Fall Meeting.
- Bhattacharyya, P., et al. (2014). **Tropical low land rice ecosystem is a net carbon sink.** *Agriculture, Ecosystems and Environment* 189: 127–135.
- Birkham, T., et al. (2014). **Near-surface water balances of waste rock dumps.** British Columbia Mine Reclamation Symposium.
- Chamberlain, S. D., et al. (2015). **Underlying ecosystem emissions exceed cattle-emitted methane from subtropical lowland pastures.** *Ecosystems* 18(6): 933–945.
- Chu, H., et al. (2014). **Net ecosystem methane and carbon dioxide exchanges in a Lake Erie coastal marsh and a nearby cropland.** *Journal of Geophysical Research: Biogeosciences* 119(5): 722–740.
- Coates, T., et al. (2014). **A year in the life: The challenges of long term methane flux measurements on a cattle-grazed landscape.** *Optical Instrumentation for Energy and Environmental Applications.* Optical Society of America.
- Emmerton, C. A., et al. (2014). **The net exchange of methane with high Arctic landscapes during the summer growing season.** *Biogeosciences* 11(12): 3095–3106.
- Fares, S. and F. Loreto (2014). **Isoprenoid emissions by the Mediterranean vegetation in castelporziano.** *Rendiconti Lincei.* 4 September 2014.
- Fleischer, E., et al. (2015). **Summer evapotranspiration in Western Siberia: a comparison between eddy covariance and Penman method formulations.** *Hydrological Processes.* 29(20): 4498–4513.
- Fortuniak, K., et al. (2014). **Singularities of the urban climate of Łódź, Central Poland.** Natural environment of Poland and its protection in Łódź University Geographical Research. Edited by E. Koboжек and T. Marszał.
- Gondwe, M. J. and W. R. Masamba (2014). **Spatial and temporal dynamics of diffusive methane emissions in the Okavango Delta, northern Botswana, Africa.** *Wetlands ecology and management* 22(1): 63–78.
- Goodrich, J., et al. (2015). **Over-riding control of methane flux temporal variability by water table dynamics in a Southern Hemisphere raised bog.** *Journal of Geophysical Research: Biogeosciences* 120: 819–831.
- Iwata, H., et al. (2014). **Cross-validation of open-path and closed-path eddy-covariance techniques for observing methane fluxes.** *Boundary-Layer Meteorology* 151(1): 95–118.
- Jha, C. S., et al. (2014). **Eddy covariance based methane flux in Sundarbans mangroves, India.** *Journal of Earth System Science* 123(5): 1089–1096.
- Knox, S. H., et al. (2015). **Agricultural peatland restoration: effects of land-use change on greenhouse gas (CO₂ and CH₄) fluxes in the Sacramento-San Joaquin Delta.** *Global Change Biology* 21(2): 750–765.
- Mammarella, I. (2015). **Eddy covariance technique: flux corrections. Workshop on EddyUH: a software for eddy covariance flux calculation.** Helsinki, 23–27 February 2015.

- Matthes, J. H., et al. (2014). **Parsing the variability in CH₄ flux at a spatially heterogeneous wetland: Integrating multiple eddy covariance towers with high-resolution flux footprint analysis.** *Journal of Geophysical Research: Biogeosciences* 119(7): 1322–1339.
- Morin, T., et al. (2014). **Environmental drivers of methane fluxes from an urban temperate wetland park.** *Journal of Geophysical Research: Biogeosciences* 119(11): 2188–2208.
- Morin, T., et al. (2014). **The seasonal and diurnal dynamics of methane flux at a created urban wetland.** *Ecological Engineering* 72: 74–83.
- Podgrajsek, E., et al. (2014). **Comparison of floating chamber and eddy covariance measurements of lake greenhouse gas fluxes.** *Biogeosciences* 11(15): 4225–4233.
- Sahlée, E., et al. (2014). **Influence from surrounding land on the turbulence measurements above a lake.** *Boundary-Layer Meteorology* 150(2): 235–258.
- Shoemaker, W., et al. (2015). **Carbon exchange between the atmosphere and subtropical forested cypress and pine wetlands.** *Biogeosciences* 12(8): 2285–2300.
- Sun, K., et al. (2014). **On-road ammonia emissions characterized by mobile, open-path measurements.** *Environmental Science and Technology* 48(7): 3943–3950.
- Tao, L., et al. (2015). **Low-power, open-path mobile sensing platform for high-resolution measurements of greenhouse gases and air pollutants.** *Applied Physics B* 119(1): 153–164.
- Torn, M. and D. B. N. Raz-Yaseef (2014). **Eddy-Covariance and auxiliary measurements, Ngee-Barrow, 2012–2013.** Next Generation Ecosystems Experiment-Arctic, Oak Ridge National Laboratory (ORNL), Oak Ridge, TN (US).
- Verkerke, J. L., et al. (2014). **Remote sensing of CO₂ leakage from geologic sequestration projects.** *International Journal of Applied Earth Observation and Geoinformation* 31: 67–77.
- Xu, L., et al. (2014). **Impact of changes in barometric pressure on landfill methane emission.** *Global Biogeochemical Cycles* 28(7): 679–695.
- Yu, Z., et al. (2014). **Dynamics of methane ebullition from a peat monolith revealed from a dynamic flux chamber system.** *Journal of Geophysical Research: Biogeosciences* 119(9): 1789–1806.
- Zhang, L. (2014). **Quantification of fugitive emissions from a biosolids lagoon.** Thesis: University of Alberta.
- Голубева, Е. and Ф. Романенко "БК 26.8 И88 ISBN No. 0-000-00000-X."
- 강남구 and 윤주열 (2015). "개회로 파장 변조 분광법과 에디 공분산 방법으로 논에서 관측된 CH₄ 플럭스 자료의 보정." *한국농림기상학회지* 17(1): 15–24.

LI-7200 Enclosed CO₂/H₂O Analyzer

- Begashaw, I., et al. (2014) **A new tool for automated data collection and complete on-site flux data processing for eddy covariance measurements.** AGU Fall Meeting.
- Birkham, T., et al. (2014). **Near-surface water balances of waste rock dumps.** British Columbia Mine Reclamation Symposium 2014.
- Buczko, U., et al. (2015). **Spatial variability at different scales and sampling requirements for in situ soil CO₂ efflux measurements on an arable soil.** *CATENA* 131: 46–55.
- Burba, G., et al. (2014). **Continuous atmospheric monitoring of the injected CO₂ behavior over geological storage sites using flux stations: latest technologies and resources.** EGU General Assembly.
- Burns, S. P., et al. (2014). **A comparison of infrared gas analyzers above a subalpine forest in complex terrain.** AGU Fall Meeting.
- Chen, Z. H., et al. (2015). **Leaf nitrogen is closely coupled to phenophases in a desert shrub ecosystem in China.** *Journal of Arid Environments* 122: 124–131.
- Cirillo, C. et al. (2015). **Artemisia: Active and Interactive Monitoring of the Forests in Protected Areas Aimed at the Sustainable Management of Nature.** EQA - International Journal of Environmental Quality 16: 35–42.
- Connan, O., et al. (2015). **In situ measurements of tritium evapotranspiration (3H-ET) flux over grass and soil using the gradient and eddy covariance experimental methods and the FAO-56 model.** *Journal of Environmental Radioactivity* 148(3): 1–9.
- De Ligne, A., et al. (2014). **IRGA gas sampling system dimensioning: Laboratory and field experiments.** 1st ICOS Conference.
- Degefe, D., et al. (2015). **Fog chemical composition and its feedback to fog water fluxes, water vapor fluxes, and microphysical evolution of two events near Paris.** *Atmospheric Research* 164–165: 328–338.

- Degeffie, D., et al. (2015). **Microphysics and energy and water fluxes of various fog types at SIRTA, France.** *Atmospheric Research* 151: 162–175.
- Drennan, W., et al. (2014). **EASI: An air–sea interaction buoy for high winds.** *Journal of Atmospheric and Oceanic Technology* 31(6): 1397–1409.
- Eder, F., et al. (2014). **Evaluation of two energy balance closure parametrizations.** *Boundary-Layer Meteorology* 151(2): 195–219.
- Eder, F., et al. (2015). **Secondary circulations at a solitary forest surrounded by semi-arid shrubland and their impact on eddy-covariance measurements.** *Agricultural and Forest Meteorology* 211–212: 115–127.
- Emmerton, C. A., et al. (2015). **Net ecosystem exchange of CO₂ with rapidly changing high Arctic landscapes.** *Global Change Biology*. First published online: 17 August 2015. DOI: 10.1111/gcb.13064.
- Emmerton, C. A., et al. (2014). **The net exchange of methane with high Arctic landscapes during the summer growing season.** *Biogeosciences* 11(12): 3095–3106.
- Fares, S. and F. Loreto (2014). **Isoprenoid emissions by the Mediterranean vegetation in Castelporziano.** *Rendiconti Lincei*. 4 September 2014.
- Fares, S. and F. Savi (2014). **Ozone dynamics in a Mediterranean Holm oak forest: comparison among transition periods characterized by different amounts of precipitation.** *Annals of Silvicultural Research* 38(1): 1–6.
- Fares, S., et al. (2014). **Simultaneous measurements of above and below canopy ozone fluxes help partitioning ozone deposition between its various sinks in a Mediterranean Oak Forest.** *Agricultural and forest meteorology* 198: 181–191.
- Fleischer, E., et al. (2015). **Summer evapotranspiration in Western Siberia: a comparison between eddy covariance and Penman method formulations.** *Hydrological Processes* 29(20): 4498–4513.
- Gentsch, L., et al. (2015). **Eddy-covariance methane flux measurements over a European beech forest.** EGU General Assembly.
- Gilmanov, T. G., et al. (2014). **Productivity and carbon dioxide exchange of leguminous crops: estimates from flux tower measurements.** *Agronomy Journal* 106(2): 545–559.
- Heiskanen, J. J., et al. (2014). **Effects of cooling and internal wave motions on gas transfer coefficients in a boreal lake.** *Tellus B* 66(1).
- Hommeltenberg, J. (2014). **Biosphere-atmosphere exchange of CO₂ and CH₄ over natural and drained bog forest ecosystems in southern Germany, München.** Dissertation: Technische Universität München.
- Hommeltenberg, J., et al. (2014). **Ecosystem scale methane fluxes in a natural temperate bog-pine forest in southern Germany.** *Agricultural and Forest Meteorology* 198: 273–284.
- Hommeltenberg, J., et al. (2014). **Can a bog drained for forestry be a stronger carbon sink than a natural bog forest?** *Biogeosciences* 11(13): 3477–3493.
- Ikawa, H., et al. (2015). **Understory CO₂, sensible heat, and latent heat fluxes in a black spruce forest in interior Alaska.** *Agricultural and Forest Meteorology* 214–215: 80–90.
- Jia, X., et al. (2014). **Biophysical controls on net ecosystem CO₂ exchange over a semiarid shrubland in northwest China.** *Biogeosciences* 11(17): 4679–4693.
- Jia, X., et al. (2015). **Energy partitioning over a semi-arid shrubland in northern China.** *Hydrological Processes*. first published online: 15 October 2015. DOI: 10.1002/hyp.10685.
- Jones, H. (2014). **Evapotranspiration, surface conductance and water-use efficiency of two young hybrid-poplar plantations in Canada's aspen parkland.** Master of science thesis: The University of British Columbia.
- Karsisto, P., et al. (2015). **Seasonal surface urban energy balance and wintertime stability simulated using three land-surface models in the high-latitude city Helsinki.** *Quarterly Journal of the Royal Meteorological Society*. First published online: 9 October 2015. DOI: 10.1002/qj.2659.
- Klemedtsson, L., et al. (2014). **Nitrous oxide emissions after sewage sludge and inorganic N-fertilization of a willow bio-energy plantation.** EGU General Assembly.
- Kulmala, L., et al. (2014). **Changes in biogeochemistry and carbon fluxes in a boreal forest after the clear-cutting and partial burning of slash.** *Agricultural and Forest Meteorology* 188: 33–44.
- Kurppa, M., et al. (2015). **Effect of seasonal variability and land use on particle number and CO₂ exchange in Helsinki, Finland.** *Urban Climate* 13: 94–109.
- Laubach, J., et al. (2015). **Combining two complementary micrometeorological methods to measure CH₄ and N₂O fluxes over pasture.** *Biogeosciences Discussions* 12(18): 15245–15299.

- Laurent, J., et al. «**FLUXALP**»: Un projet de développement d'une station de mesures éco-climatiques au Col Du Lautaret, Hautes-Alpes, France.
- Lohila, A., et al. (2015). **Carbon dioxide and energy fluxes over a northern boreal lake.** *Boreal Environment Research* 20(4): 474–488.
- Loogus, M. (2014). **Süsinikuvoo mõõtmised turbulentse kovariatsiooni meetodil.**
- Lund, M., et al. (2015). **Low impact of dry conditions on the CO₂ exchange of a Northern-Norwegian blanket bog.** *Environmental Research Letters* 10(2): 025004.
- Mammarella, I., et al. (2015). **Carbon dioxide and energy fluxes over a small boreal lake in Southern Finland.** *Journal of Geophysical Research: Biogeosciences* 120(7): 1296–1314.
- Mammarella, I. (2015). **Eddy Covariance technique: flux corrections.** Workshop on EddyUH: a software for eddy covariance flux calculation. Helsinki.
- Marcolla, B., et al. (2014). **Methods and uncertainties in the experimental assessment of horizontal advection.** *Agricultural and Forest Meteorology* 198: 62–71.
- Mauder, M., et al. (2013). **A strategy for quality and uncertainty assessment of long-term eddy-covariance measurements.** *Agricultural and Forest Meteorology* 169: 122–135.
- Meshalkina, J., et al. (2015). **Central Russia agroecosystem monitoring with CO₂ fluxes analysis by eddy covariance method.** Abstract book, 9th International Soil Science Congress.
- Moat, B., et al. (2015). **Metadata for the WAGES instrumentation deployed on the RRS James Clark Ross between May 2010 and September 2013.** Southampton, National Oceanography Centre.
- Nordbo, A., et al. (2014). **Sorption-Caused Attenuation and Delay of Water Vapor Signals in Eddy-Covariance Sampling Tubes and Filters.** *Journal of Atmospheric and Oceanic Technology* 31(12): 2629–2649.
- Nordbo, A., et al. (2015). **Urban surface cover determined with airborne lidar at 2m resolution - Implications for surface energy balance modelling.** *Urban Climate* 13: 52–72.
- Oechel, W. C., et al. (2014). **Annual patterns and budget of CO₂ flux in an Arctic tussock tundra ecosystem.** *Journal of Geophysical Research: Biogeosciences* 119(3): 323–339.
- Ortiz, P. S., et al. (2015). **CO₂ and CH₄ exchange by Phragmites australis under different climates.** EGU General Assembly.
- O'Shea, S., et al. (2014). **Methane and carbon dioxide fluxes and their regional scalability for the European Arctic wetlands during the MAMM project in summer 2012.** *Atmospheric Chemistry and Physics* 14(23): 13159–13174.
- Pelletier, L., et al. (2015). **Can boreal peatlands with pools be net sinks for CO₂?** *Environmental Research Letters* 10(3): 035002.
- Pelletier, L., et al. (2014). **Effect of open water pools on ecosystem scale surface-atmosphere carbon dioxide exchange in a boreal peatland.** *Biogeochemistry* 124(1–3): 291–304.
- Shoemaker, J. K., et al. (2014). **Forest ecosystem changes from annual methane source to sink depending on late summer water balance.** *Geophysical Research Letters* 41(2): 673–679.
- Sievers, J., et al. (2014). **Estimating local atmosphere-surface fluxes using eddy covariance and numerical Ogive optimization.** *Atmospheric Chemistry and Physics Discussions* 14(15): 21387–21432.
- Sievers, J., et al. (2015). **Estimating surface fluxes using eddy covariance and numerical ogive optimization.** *Atmospheric Chemistry and Physics* 15(4): 2081–2103.
- Sievers, J., et al. (2015). **Winter observations of CO₂ exchange between sea-ice and the atmosphere in a coastal fjord environment.** *The Cryosphere Discussions* 9(1): 45–75.
- Unit, T. A. (2014). **Two bogs in the Canadian Hudson Bay lowlands and a temperate bog reveal similar annual net ecosystem exchange of CO₂.** *Arctic, Antarctic, and Alpine Research* 46(1): 103–113.
- Welch, C. M., et al. (2015). **The impacts of mountain pine beetle disturbance on the energy balance of snow during the melt period.** *Hydrological Processes*. First published online: 13 September 2015. DOI: 10.1002/hyp.10638.
- Winderlich, J., et al. (2014). **Inferences from CO₂ and CH₄ concentration profiles at the Zotino Tall Tower Observatory (ZOTTO) on regional summertime ecosystem fluxes.** *Biogeosciences* 11(7): 2055–2068.
- Xie, J., et al. (2015). **Irregular precipitation events in control of seasonal variations in CO₂ exchange in a cold desert-shrub ecosystem in northwest China.** *Journal of Arid Environments* 120: 33–41.

LI-7500/A Open Path CO₂/H₂O Analyzer

- Abraha, M., et al. (2014). **Evapotranspiration of annual and perennial biofuel crops in a variable climate.** *GCB Bioenergy*.
- Aguilos, M., et al. (2014). **Dynamics of ecosystem carbon balance recovering from a clear-cutting in a cool-temperate forest.** *Agricultural and Forest Meteorology* 197: 26–39.
- Alberto, M. C. R., et al. (2014). **Actual evapotranspiration and dual crop coefficients for dry-seeded rice and hybrid maize grown with overhead sprinkler irrigation.** *Agricultural Water Management* 136: 1–12.
- Alberto, M. C. R., et al. (2014). **Measuring methane flux from irrigated rice fields by eddy covariance method using open-path gas analyzer.** *Field Crops Research* 160: 12–21.
- Amatya, P. M., et al. (2015). **Mapping regional distribution of land surface heat fluxes on the southern side of the central Himalayas using TESEBS.** *Theoretical and Applied Climatology* 23 April 2015. DOI: 10.1007/s00704-015-1466-2.
- Andersson, A., et al. (2014). **Using a high-frequency fluorescent oxygen probe in atmospheric eddy covariance applications.** *Journal of Atmospheric and Oceanic Technology* 31(11): 2498–2511.
- Artigas, F., et al. (2015). **Long term carbon storage potential and CO₂ sink strength of a restored salt marsh in New Jersey.** *Agricultural and Forest Meteorology* 200: 313–321.
- Babel, W., et al. (2014). **Pasture degradation modifies the water and carbon cycles of the Tibetan highlands.** *Biogeosciences Discussions* 11(6): 8861–8923.
- Bai, J., et al. (2015). **Seasonal and inter-annual variations in carbon fluxes and evapotranspiration over cotton field under drip irrigation with plastic mulch in an arid region of Northwest China.** *Journal of Arid Land* 7(2): 272–284.
- Balogh, J., et al. (2015). **Soil CO₂ efflux and production rates as influenced by evapotranspiration in a dry grassland.** *Plant and Soil* 388: 157–173.
- Bao, X., et al. (2014). **Interannual variation in carbon sequestration depends mainly on the carbon uptake period in two croplands on the North China Plain.** *PLoS one* 9(10): e110021.
- Barr, J. G., et al. (2014). **Seasonal evapotranspiration patterns in mangrove forests.** *Journal of Geophysical Research: Atmospheres* 119(7): 3886–3899.
- Bawazir, A. S., et al. (2014). **Assessment of the crop coefficient for saltgrass under native riparian field conditions in the desert southwest.** *Hydrological Processes* 28(25): 6163–6171.
- Begashaw, I., et al. (2014) **A new tool for automated data collection and complete on-site flux data processing for eddy covariance measurements.** AGU Fall Meeting.
- Bhattacharyya, P., et al. (2014). **Tropical low land rice ecosystem is a net carbon sink.** *Agriculture, Ecosystems and Environment* 189: 127–135.
- Biederman, J., et al. (2014). **Increased evaporation following widespread tree mortality limits streamflow response.** *Water Resources Research* 50(7): 5395–5409.
- Billesbach, D., et al. (2014). **Growing season eddy covariance measurements of carbonyl sulfide and CO₂ fluxes: COS and CO₂ relationships in Southern Great Plains winter wheat.** *Agricultural and Forest Meteorology* 184: 48–55.
- Biudes, M. S., et al. (2014). **Modelling gross primary production of a tropical semi-deciduous forest in the southern Amazon Basin.** *International Journal of Remote Sensing* 35(4): 1540–1562.
- Bohnenstengel, S., et al. (2015). **Meteorology, air quality, and health in London: The ClearLo project.** *Bulletin of the American Meteorological Society*. May 2015.
- Boylan, P., et al. (2014). **Characterization and mitigation of water vapor effects in the measurement of ozone by chemiluminescence with nitric oxide.** *Atmospheric Measurement Techniques* 7(5): 1231–1244.
- Brown, M. G., et al. (2014). **Evapotranspiration and canopy characteristics of two lodgepole pine stands following mountain pine beetle attack.** *Hydrological Processes* 28(8): 3326–3340.
- Brown, S., et al. (2014). **Atmospheric and soil moisture controls on evapotranspiration from above and within a Western Boreal Plain aspen forest.** *Hydrological Processes* 28(15): 4449–4462.
- Budishchev, A., et al. (2014). **Evaluation of a plot-scale methane emission model using eddy covariance observations and footprint modelling.** *Biogeosciences* 11(17): 4651–4664.

- Burba, G., et al. (2014). **Advancements in micro-meteorological technique for monitoring CH₄ release from remote permafrost regions: Principles, emerging research, and latest updates.** EGU General Assembly.
- Burba, G., et al. (2014). **Continuous atmospheric monitoring of the injected CO₂ behavior over geological storage sites using flux stations: latest technologies and resources.** EGU General Assembly.
- Burns, S. P., et al. (2014) **A comparison of infrared gas analyzers above a subalpine forest in complex terrain.** AGU Fall Meeting.
- Byun, K., et al. (2014). **Dual-model approaches for evapotranspiration analyses over homo-and heterogeneous land surface conditions.** *Agricultural and Forest Meteorology* 197: 169–187.
- Cammalleri, C., et al. (2014). **Mapping daily evapotranspiration at field scales over rainfed and irrigated agricultural areas using remote sensing data fusion.** *Agricultural and Forest Meteorology* 186: 1–11.
- Campbell, D. I., et al. (2015). **Variations in CO₂ exchange for dairy farms with year-round rotational grazing on drained peatlands.** *Agriculture, Ecosystems and Environment* 202: 68–78.
- Carrasco-Benavides, M., et al. (2014). **Parameterization of the satellite-based model (METRIC) for the estimation of instantaneous surface energy balance components over a drip-irrigated vineyard.** *Remote Sensing* 6(11): 11342–11371.
- Chamberlain, S. D., et al. (2015). **Underlying ecosystem emissions exceed cattle-emitted methane from subtropical lowland pastures.** *Ecosystems* 18(6): 933–945.
- Chang, K.H., et al. (2014). **A simple crop phenology algorithm in the land surface model CN-CLASS.** *Agronomy Journal* 106(1): 297–308.
- Charuchittipan, D., et al. (2014). **Extension of the averaging time in eddy-covariance measurements and its effect on the energy balance closure.** *Boundary-Layer Meteorology* 152 (3): 303–327.
- Chen, F., et al. (2015). **An observational and modeling study of impacts of bark beetle-caused tree mortality on surface energy and hydrological cycles.** *Journal of Hydrometeorology* 16(2): 744–761.
- Chen, J., et al. (2015). **Ergodicity test of the eddy-covariance technique.** *Atmospheric Chemistry and Physics* 15(17): 9929–9944.
- Chow, W. T., et al. (2014). **Seasonal dynamics of a suburban energy balance in Phoenix, Arizona.** *International Journal of Climatology* 34(15): 3863–3880.
- Chowdhuri, S., et al. (2015). **Relationship between the momentum and scalar fluxes close to the ground during the indian post-monsoon period.** *Boundary-Layer Meteorology* 154(2): 333–348.
- Christen, A., et al. (2014). **Quantifying summertime greenhouse gases fluxes from soils in various stages of restoration in the Burns Bog Ecological Conservancy Area.** University of British Columbia. DOI: 10.13140/2.1.3977.5042.
- Churakova, O. V., et al. (2014). **Increasing relevance of spring temperatures for Norway spruce trees in Davos, Switzerland, after the 1950s.** *Trees* 28(1): 183–191.
- Coates, T., et al. (2014). **A year In the life: The challenges of long term methane flux measurements on a cattle-grazed landscape.** Optical Instrumentation for Energy and Environmental Applications. Optical Society of America.
- Crawford, B. and A. Christen (2015). **Spatial source attribution of measured urban eddy covariance CO₂ fluxes.** *Theoretical and Applied Climatology* 119 (3–4): 733–755.
- Crawford, B. and A. Christen (2014). **Spatial variability of carbon dioxide in the urban canopy layer and implications for flux measurements.** *Atmospheric Environment* 98: 308–322.
- Crilley, L., et al. (2014). **Sources and contributions of wood smoke during winter in London: assessing local and regional influences.** *Atmospheric Chemistry and Physics Discussions* 14 (19): 27459–27530.
- Crilley, L., et al. (2015). **Sources and contributions of wood smoke during winter in London: assessing local and regional influences.** *Atmospheric Chemistry and Physics* 15(6): 3149–3171.
- Dadi, T., et al. (2015). **Post-wildfire effects on carbon and water vapour dynamics in a Spanish black pine forest.** *Environmental Science and Pollution Research* 22(7): 4851–4862.
- Degefie, D., et al. (2015). **Microphysics and energy and water fluxes of various fog types at SIRT, France.** *Atmospheric Research* 151: 162–175.

- Della Chiesa, S., et al. (2014). **Modelling changes in grassland hydrological cycling along an elevational gradient in the Alps.** *Ecohydrology* 7 (6): 1453–1473.
- DeLucia, N., et al. (2015). **Management practices and controls on methane emissions from subtropical wetlands.** EGU General Assembly.
- Diaz, M. and D. Roberti (2015). **Calculation of carbon exchanges in rice agroecosystem with application of Gap-filling techniques.** *Ciência e Natura*.
- Diaz, M. and D. Roberti (2015). **Cálculo das trocas de carbono num agroecossistema de arroz com aplicação de técnicas de preenchimento de falhas.** *Ciência e Natura*.
- Drennan, W., et al. (2014). **EASI: An air–sea interaction buoy for high winds.** *Journal of Atmospheric and Oceanic Technology* 31(6): 1397–1409.
- Du, Q., et al. (2014). **Effects of different gap filling methods and land surface energy balance closure on annual net ecosystem exchange in a semiarid area of China.** *Science China Earth Sciences* 57(6): 1340–1351.
- Eder, F., et al. (2014). **Evaluation of two energy balance closure parametrizations.** *Boundary-Layer Meteorology* 151(2): 195–219.
- Eder, F., et al. (2015). **Mesoscale eddies affect near-surface turbulent exchange: evidence from lidar and tower measurements.** *Journal of Applied Meteorology and Climatology* 54(1): 189–206.
- El-Madany, T., et al. (2014). **Low-level jets and above-canopy drainage as causes of turbulent exchange in the nocturnal boundary layer.** *Biogeosciences* 11(16): 4507–4519.
- Else, B., et al. (2014). **Surface energy budget of landfast sea ice during the transitions from winter to snowmelt and melt pond onset: The importance of net longwave radiation and cyclone forcings.** *Journal of Geophysical Research: Oceans* 119(6): 3679–3693.
- Emmel, C. (2014). **Vertical distribution of carbon dioxide, water vapour, momentum and energy exchange within and above a forest stand affected by the mountain pine beetle.** Doctoral dissertation: University of British Columbia. Vancouver, Canada.
- Emmel, C., et al. (2014). **Vertical distribution of carbon dioxide sources and sinks in a recovering mountain pine beetle-attacked lodgepole pine stand.** *Agricultural and Forest Meteorology* 195: 108–122.
- Emmerton, C. A., et al. (2014). **The net exchange of methane with high Arctic landscapes during the summer growing season.** *Biogeosciences* 11(12): 3095–3106.
- Etheridge, D., et al. (2014) **Metadata report: Arcturus atmospheric greenhouse gas monitoring.** Geoscience Australia and CSIRO.
- Evrendilek, F. (2014). **Assessing CO₂ sink/source strength of a degraded temperate peatland: atmospheric and hydrological drivers and responses to extreme events.** *Ecohydrology*. First published online: 22 December 2014. DOI: 10.1002/eco.1592.
- Evrendilek, F. (2014). **Assessing neural networks with wavelet denoising and regression models in predicting diel dynamics of eddy covariance-measured latent and sensible heat fluxes and evapotranspiration.** *Neural Computing and Applications* 24(2): 327–337.
- Evrendilek, F. (2014). **Modeling net ecosystem carbon dioxide exchange using temporal neural networks after wavelet denoising.** *Geographical Analysis* 46(1): 37–52.
- Fares, S. and F. Loreto (2014). **Isoprenoid emissions by the Mediterranean vegetation in castelporziano.** *Rendiconti Lincei*. 4 September 2014.
- Flerchinger, G. and M. Seyfried (2014). **Comparison of methods for estimating evapotranspiration in a small rangeland catchment.** *Vadose Zone Journal* 13(4).
- Font, A., et al. (2015). **Daytime CO₂ urban surface fluxes from airborne measurements, eddy-covariance observations and emissions inventory in Greater London.** *Environmental Pollution* 196: 98–106.
- Frank, J. M., et al. (2014). **Ecosystem CO₂/H₂O fluxes are explained by hydraulically limited gas exchange during tree mortality from spruce bark beetles.** *Journal of Geophysical Research: Biogeosciences* 119(6): 1195–1215.
- Fratini, G. and M. Mauder (2014). **Towards a consistent eddy-covariance processing: an intercomparison of EddyPro and TK3.** *Atmospheric Measurement Techniques* 7(7): 2273–2281.

- Gao, S., et al. (2015). **Ecosystem carbon (CO₂ and CH₄) fluxes of a *Populus deltoides* plantation in subtropical China during and post clear-cutting.** *Forest Ecology and Management* 357: 206–219.
- Gao, Y., et al. (2015). **Evapotranspiration over artificially planted shrub communities in the shifting sand dune area of the Tengger Desert, north central China.** *Ecohydrology*. First published online: 12 May 2015.
- Gaofeng, Z., et al. (2014). **Energy flux partitioning and evapotranspiration in a sub-alpine spruce forest ecosystem.** *Hydrological Processes* 28(19): 5093–5104.
- Garcia, M., et al. (2014). **Accuracy of the temperature-vegetation dryness Index using MODIS under water-limited vs. energy-limited evapotranspiration conditions.** *Remote Sensing of Environment* 149: 100–117.
- Gilmanov, T. G., et al. (2014). **Productivity and carbon dioxide exchange of leguminous crops: estimates from flux tower measurements.** *Agronomy Journal* 106(2): 545–559.
- Glenn, E. P., et al. (2015). **Wide-area ratios of evapotranspiration to precipitation in monsoon-dependent semiarid vegetation communities.** *Journal of Arid Environments* 117: 84–95.
- Goldbach, A. and W. Kuttler (2014). **Turbulent heat fluxes above a suburban reservoir—a case study from Germany.** *Journal of Hydrometeorology* 16: 244–260.
- Gong, J., et al. (2014). **Climatic sensitivity of the CO₂ flux in a cutaway boreal peatland cultivated with a perennial bioenergy crop (*Phalaris arundinaceae*, L.): Beyond diplotelmic modeling.** *Agricultural and Forest Meteorology* 198: 232–249.
- González, A. Z. (2014). **Efecto del sistema de riego y clima en la eficiencia del uso de agua de nogal pecanero.** *REVISTA TERRA LATINOAMERICANA* 32 (1): 23–34.
- Goodrich, J., et al. (2015). **Overriding control of methane flux temporal variability by water table dynamics in a Southern Hemisphere raised bog.** *Journal of Geophysical Research: Biogeosciences* 120(5): 819–831.
- Gou, Q., et al. (2015) **Microclimate and CO₂ fluxes on continuous fine days in the Xihu desert wetland, China.** *Journal of Arid Land*: 7(3): 318–327.
- Gou, Q., et al. (2014). **Analysis of the microclimate and CO₂ flux characteristics in arid desert wetland ecosystem in summer: A case study in Xihu desert wetland, Dunhuang, China.** *Journal of Earth System Science* 123(7): 1491–1499.
- Graf, A., et al. (2014). **Spatiotemporal relations between water budget components and soil water content in a forested tributary catchment.** *Water Resources Research* 50(6): 4837–4857.
- Gu, L., et al. (2015). **Comparison of the surface energy budget between regions of seasonally frozen ground and permafrost on the Tibetan Plateau.** *Atmospheric Research* 153: 553–564.
- Guo, W., et al. (2014). **Variation of NEE and its affecting factors in a vineyard of arid region of northwest China.** *Atmospheric Environment* 84: 349–354.
- Guo, X. F., et al. (2014). **Urban boundary-layer stability and turbulent exchange during consecutive episodes of particle air pollution in Beijing, China.** *Atmospheric and Oceanic Science Letters* 7: 62–66.
- Hamerlynck, E. P., et al. (2014). **Water use efficiency of annual-dominated and bunchgrass-dominated savanna intercanopy space.** *Ecohydrology* 7(4): 1208–1215.
- Han, C., et al. (2014). **Estimates of effective aerodynamic roughness length over mountainous areas of the Tibetan Plateau.** *Quarterly Journal of the Royal Meteorological Society* 141(689): 1457–1465.
- Han, G., et al. (2014). **Agricultural reclamation effects on ecosystem CO₂ exchange of a coastal wetland in the Yellow River Delta.** *Agriculture, Ecosystems and Environment* 196: 187–198.
- Han, G., et al. (2014). **Ecosystem photosynthesis regulates soil respiration on a diurnal scale with a short-term time lag in a coastal wetland.** *Soil Biology and Biochemistry* 68: 85–94.
- Hanis, K. L., et al. (2015). **Carbon exchange over four growing seasons for a subarctic sedge fen in northern Manitoba, Canada.** *Arctic Science*. Published online: 21 September 2015.
- Hauwert, N. M. and J. M. Sharp (2014). **Measuring autogenic recharge over a karst aquifer utilizing eddy covariance evapotranspiration.** *Journal of Water Resource and Protection* 6: 869–879.
- Heilman, J. L., et al. (2014). **Water storage capacity controls energy partitioning and water use in karst ecosystems on the Edwards Plateau, Texas.** *Ecohydrology* 7(1): 127–138.

- Hilker, T., et al. (2014). **Potentials and limitations for estimating daytime ecosystem respiration by combining tower-based remote sensing and carbon flux measurements.** *Remote Sensing of Environment* 150: 44–52.
- Hiller, R. V., et al. (2014). **Aircraft-based CH₄ flux estimates for validation of emissions from an agriculturally dominated area in Switzerland.** *Journal of Geophysical Research: Atmospheres* 119 (8): 4874–4887.
- Hirano, T., et al. (2015). **Evapotranspiration of tropical peat swamp forests.** *Global Change Biology* 21 (5):1914–1927.
- Hommeltenberg, J. (2014). **Biosphere-atmosphere exchange of CO₂ and CH₄ over natural and drained bog forest ecosystems in southern Germany.** Dissertation: München, Technische Universität München.
- Hong, J., et al. (2014). **Wet surface resistance of forest canopy in monsoon Asia: Implications for eddy-covariance measurement of evapotranspiration.** *Hydrological Processes* 28(1): 37–42.
- Hoppe, C., et al. (2014). **A variational data assimilation system for soil–atmosphere flux estimates for the Community Land Model (CLM3. 5).** *Geoscientific Model Development* 7(3): 1025–1036.
- Hoppmann, M., et al. (2015). **Ice platelets below Weddell Sea landfast sea ice.** *Annals of Glaciology* 56(69): 175.
- Hu, Z., et al. (2014). **Partitioning of evapotranspiration through oxygen isotopic measurements of water pools and fluxes in a temperate grassland.** *Journal of Geophysical Research: Biogeosciences* 119(3): 358–372.
- Huang, H., et al. (2015). **Atmospheric boundary layer structure and turbulence during sea fog on the southern China coast.** *Monthly Weather Review* 143: 1907–1923.
- Huang, L. and X. Wen (2014). **Temporal variations of atmospheric water vapor δD and $\delta^{18}O$ above an arid artificial oasis cropland in the Heihe River Basin.** *Journal of Geophysical Research: Atmospheres* 119(19): 11,456–11,476.
- Hughes, D., et al. (2014). **Using targeted short term field investigations to calibrate and evaluate the structure of a hydrological model.** *Hydrological Processes* 28(5): 2794–2809.
- Hunt, E. D., et al. (2014). **Monitoring the effects of rapid onset of drought on non-irrigated maize with agronomic data and climate-based drought indices.** *Agricultural and Forest Meteorology* 191: 1–11.
- Huo, Q., et al. (2014). **Inference of emission rate using the inverse-dispersion method for the multi-source problem.** *Agricultural and Forest Meteorology* 191: 12–21.
- Huo, Q., et al. (2015). **Effects of surface source/sink distributions on the flux–gradient similarity in the unstable surface layer.** *Theoretical and Applied Climatology* 119(1–2): 313–322.
- Huo, Q., et al. (2015). **Estimating ammonia emissions from a winter wheat cropland in North China Plain with field experiments and inverse dispersion modeling.** *Atmospheric Environment* 104: 1–10.
- Igarashi, Y., et al. (2015). **Environmental control of canopy stomatal conductance in a tropical deciduous forest in northern Thailand.** *Agricultural and Forest Meteorology* 202: 1–10.
- Ikawa, H. and W. C. Oechel (2014). **Spatial and temporal variability of air-sea CO₂ exchange of alongshore waters in summer near Barrow, Alaska.** *Estuarine, Coastal and Shelf Science* 141: 37–46.
- Ikawa, H. and W. C. Oechel (2015). **Temporal variations in air sea CO₂ exchange near large kelp beds near San Diego, California.** *Journal of Geophysical Research: Oceans* 120(1): 50–63.
- Imukova, K., et al. (2015). **Energy balance closure on a winter wheat stand: Comparing the eddy covariance technique with the soil water balance method.** *Biogeosciences Discussions* 12(9): 6783–6820.
- Isabelle, P., et al. (2015). **Applicability of the bulk-transfer approach to estimate evapotranspiration from boreal peatlands.** *Journal of Hydrometeorology*. 16(4): 1521–1539.
- Iwata, H., et al. (2014). **Cross-validation of open-path and closed-path eddy-covariance techniques for observing methane fluxes.** *Boundary-Layer Meteorology* 151(1): 95–118.
- Jacobs, C., et al. (2015). **Assessment of evaporative water loss from Dutch cities.** *Building and Environment* 83: 27–38.
- Järvi, L., et al. (2014). **Development of the surface urban energy and water balance scheme (SUEWS) for cold climate cities.** *Geoscientific Model Development* 7(4): 1691–1711.

- Jiang, G., et al. (2014). **Analysis of light use efficiency and gross primary productivity based on remote sensing data over a phragmites-dominated wetland in Zhangye, China.** SPIE Asia Pacific Remote Sensing, International Society for Optics and Photonics.
- Jiang, Y., et al. (2014). **Dynamics of carbon fluxes with responses to vegetation, meteorological and terrain factors in the south-eastern Tibetan Plateau.** *Environmental Earth Sciences* 72(11): 4551–4565.
- Jing, C. Q., et al. (2014). **Comparison of root water uptake functions to simulate surface energy fluxes within a deep-rooted desert shrub ecosystem.** *Hydrological Processes* 28(21): 5436–5449.
- Jing, Y., et al. (2014). **Carbon dioxide fluxes over a temperate meadow in eastern Inner Mongolia, China.** *Environmental Earth Sciences* 72(11): 4401–4411.
- Jones, H. (2014). **Evapotranspiration, surface conductance and water-use efficiency of two young hybrid-poplar plantations in Canada's aspen parkland.** Thesis: The University of British Columbia.
- Kalogridis, C., et al. (2014). **Concentrations and fluxes of isoprene and oxygenated VOCs at a French Mediterranean oak forest.** *Atmospheric Chemistry and Physics* 14(18): 10085–10102.
- Kang, M., et al. (2015). **Responses of energy partitioning and surface resistance to drought in a poplar plantation in northern China.** *Biogeosciences Discussions* 12(1): 345–388.
- Kang, X., et al. (2014). **Modeling carbon fluxes using multi-temporal MODIS Imagery and CO₂ eddy flux tower data in Zoige Alpine Wetland, South-West China.** *Wetlands* 34(3): 603–618.
- Katata, G., et al. (2014). **A numerical study of the effects of aerosol hygroscopic properties to dry deposition on a broad-leaved forest.** *Atmospheric Environment* 97: 501–510.
- Khomik, M., et al. (2014). **On the causes of rising gross ecosystem productivity in a regenerating clearcut environment: leaf area vs. species composition.** *Tree Physiology* 34(7): 686–700.
- Kirschbaum, M. U., et al. (2015). **Modelling carbon and water exchange of a grazed pasture in New Zealand constrained by eddy covariance measurements.** *Science of The Total Environment* 512: 273–286.
- Knox, S. H., et al. (2014). **Agricultural peatland restoration: effects of land-use change on greenhouse gas (CO₂ and CH₄) fluxes in the Sacramento-San Joaquin Delta.** *Global Change Biology* 21(2): 750–765.
- Kongoli, C., et al. (2014). **Evaluation of a two-source snow-vegetation energy balance model for estimating surface energy fluxes in a rangeland ecosystem.** *Journal of Hydrometeorology* 15(1): 143–158.
- Kool, D., et al. (2014). **Spatial and diurnal below canopy evaporation in a desert vineyard: Measurements and modeling.** *Water Resources Research* 50(8): 7035–7049.
- Kotani, A., et al. (2014). **Temporal variations in the linkage between the net ecosystem exchange of water vapour and CO₂ over boreal forests in eastern Siberia.** *Ecohydrology* 7(2): 209–225.
- Krauss, K. W., et al. (2015). **Approximations of stand water use versus evapotranspiration from three mangrove forests in southwest Florida, USA.** *Agricultural and Forest Meteorology* 213: 291–303.
- Krofcheck, D. (2015). **Bridging structure and function in semi-arid ecosystems by integrating remote sensing and ground based measurements.** Dissertation: University of New Mexico.
- Krofcheck, D. J., et al. (2014). **Detecting mortality induced structural and functional changes in a piñon-juniper woodland using Landsat and RapidEye time series.** *Remote Sensing of Environment* 151: 102–113.
- Kulmala, L., et al. (2014). **Changes in biogeochemistry and carbon fluxes in a boreal forest after the clear-cutting and partial burning of slash.** *Agricultural and Forest Meteorology* 188: 33–44.
- Lagomasino, D., et al. (2015). **Spatial and temporal variability in spectral-based surface energy evapotranspiration measured from Landsat 5TM across two mangrove ecotones.** *Agricultural and Forest Meteorology*. 213: 304–316.
- Lapina, L. E., et al. (2015). **The Factor Analysis of Meso-Oligotrophic Peatland Medla-Pev-Njur Data.** *Procedia Earth and Planetary Science* 15(9): 688–693.
- Laurent, J., et al. «FLUXALP»: **Un projet de développement d'une station de mesures éco-climatiques au Col Du Lautaret, Hautes-Alpes, France.**

- Lee, S.C., et al. (2015). **Investigating effect of environmental controls on dynamics of CO₂ budget in a subtropical estuarial marsh wetland ecosystem.** *Environmental Research Letters* 10(2): 025005.
- Lee, T. R., et al. (2015). **Meteorological controls on the diurnal variability of carbon monoxide mixing ratio at a mountaintop monitoring site in the Appalachian Mountains.** *Tellus B*: 67(1).
- Levy, P. E., and A. Gray (2015). **Greenhouse gas balance of a semi-natural peatbog in northern Scotland.** *Environmental Research Letters* 10(9): 94019.
- Lewicki, J. and G. Hilley (2014). **Multi-scale observations of the variability of magmatic CO₂ emissions, Mammoth Mountain, CA, USA.** *Journal of Volcanology and Geothermal Research* 284: 1–15.
- Li, D., et al. (2015). **Contrasting responses of urban and rural surface energy budgets to heat waves explain synergies between urban heat islands and heat waves.** *Environmental Research Letters* 10 (5).
- Li, D., et al. (2015). **The k⁻¹ scaling of air temperature spectra in atmospheric surface layer flows.** *Quarterly Journal of the Royal Meteorological Society*. Published Online: 16 September 2015. DOI: 10.1002/qj.2668.
- Li, F. and B. Wang (2015). **A study of fitting a swamp meadow ecosystem evapotranspiration to a model based on the Penman-Monteith equation.** *Journal of Chemistry* 2015(1): 1–8.
- Li, H., et al. (2014). **Seasonal and interannual variations of ecosystem photosynthetic features in an alpine dwarf shrubland on the Qinghai-Tibetan Plateau, China.** *Photosynthetica* 52(3): 321–331.
- Li, L., et al. (2014). **Growing season net ecosystem CO₂ exchange of two desert ecosystems with alkaline soils in Kazakhstan.** *Ecology and Evolution* 4(1): 14–26.
- Li, M., et al. (2014). **A 3-year dataset of sensible and latent heat fluxes from the Tibetan Plateau, derived using eddy covariance measurements.** *Theoretical and Applied Climatology*. 28 October 2014.
- Li, N., and Jiang, L. (2015). **Temporal variations in the surface soil heat flux over maize and grass surfaces in northwest China.** *Soil Science Society of America Journal* 79(4): 1084–1093.
- Li, X., et al. (2014). **Ecological restoration and recovery in the wind-blown sand hazard areas of northern China: relationship between soil water and carrying capacity for vegetation in the Tengger Desert.** *Science China Life Sciences* 57(5): 539–548.
- Li, Z., et al. (2015). **Turbulent transfer coefficient and roughness length in a high-altitude lake, Tibetan Plateau.** *Theoretical and Applied Climatology*. 17 April 2015.
- Lietzke, B., et al. (2015). **On the controlling factors for the variability of carbon dioxide flux in a heterogeneous urban environment.** *International Journal of Climatology*. First published online: 24 January 2015.
- Lindauer, M., et al. (2014). **Net ecosystem exchange over a non-cleared wind-throw-disturbed upland spruce forest—Measurements and simulations.** *Agricultural and Forest Meteorology* 197: 219–234.
- Liu, C., et al. (2015). **Optimization and evaluation of vegetation photosynthesis and respiration model using the measurements collected from the forest site of subtropical coniferous-evergreen.** *Chinese Journal of Plant Ecology* 39: 388–397.
- Liu, H., et al. (2015). **Eddy covariance measurements of water vapor and CO₂ fluxes above the Erhai Lake.** *Science China Earth Sciences* 58(3): 317–328.
- Liu, X., et al. (2015). **Partitioning evapotranspiration in an intact forested watershed in southern China.** *Ecohydrology*. 8(6): 1037–1047.
- Logsdon, S. D., et al. (2014). **Comparison of corn transpiration, eddy covariance, and soil water loss.** *Soil Science Society of America Journal* 78(4): 1214–1223.
- Lund, M., et al. (2015). **Low impact of dry conditions on the CO₂ exchange of a Northern-Norwegian blanket bog.** *Environmental Research Letters* 10(2): 25004.
- Lv, L. (2014). **Linking montane soil moisture measurements to evapotranspiration using inverse numerical modeling.** Dissertation: Utah State University. Paper 3323.
- Ma, N., et al. (2014). **Observation of mega-dune evaporation after various rain events in the hinterland of Badain Jaran Desert, China.** *Chinese Science Bulletin* 59(2): 162–170.
- Ma, Y., et al. (2015). **Atmospheric boundary layer characteristics and land-atmosphere energy transfer in the Third Pole area.** *Iahs-aish Proceedings and Reports* 368: 27–32.

- Malone, S. L. (2014). **Hydrology drives everglades ecosystem function: implications for ecosystem vulnerability to drought, energy balance, climate teleconnections and climate change.** Dissertation, The University of Alabama Tuscaloosa.
- Malone, S. L., et al. (2014). **El Niño Southern Oscillation (ENSO) enhances CO₂ exchange rates in freshwater marsh ecosystems in the Florida Everglades.** *PLoS one* 9(12): e115058.
- Malone, S. L., et al. (2014). **Seasonal patterns in energy partitioning of two freshwater marsh ecosystems in the Florida Everglades.** *Journal of Geophysical Research: Biogeosciences* 119(8): 1487–1505.
- Mamadou, O., et al. (2014). **Energy fluxes and surface characteristics over a cultivated area in Benin: daily and seasonal dynamics.** *Hydrology and Earth System Sciences* 18(3): 893–914.
- Mammarella, I. (2015). **Eddy Covariance technique: flux corrections.** Workshop on EddyUH: a software for eddy covariance flux calculation. Helsinki, 23–27 February 2015.
- Marcolla, B., et al. (2014). **Methods and uncertainties in the experimental assessment of horizontal advection.** *Agricultural and Forest Meteorology* 198: 62–71.
- Masseroni, D., et al. (2014). **Limitations and improvements of the energy balance closure with reference to experimental data measured over a maize field.** *Atmosfera* 27(4): 335–352.
- Matsuura, S., et al. (2014). **Seasonal carbon dynamics and the effects of manure application on carbon budget of a managed grassland in a temperate, humid region in Japan.** *Grassland Science* 60(2): 76–91.
- Matteucci, M., et al. (2014). **Soil and ecosystem carbon fluxes in a Mediterranean forest during and after drought.** *Agrochimica* 58(1): 91–115.
- Matthes, J. H., et al. (2014). **Parsing the variability in CH₄ flux at a spatially heterogeneous wetland: Integrating multiple eddy covariance towers with high-resolution flux footprint analysis.** *Journal of Geophysical Research: Biogeosciences* 119(7): 1322–1339.
- Matthes, J., et al. (2015). **Predicting landscape-scale CO₂ flux at a pasture and rice paddy with long-term hyperspectral canopy reflectance measurements.** *Biogeosciences Discussion* 12: 5079–5122.
- Mauder, M., et al. (2013) **A strategy for quality and uncertainty assessment of long-term eddy-covariance measurements.** *Agricultural and Forest Meteorology* 169: 122–135
- Mbufong, H., et al. (2014). **Assessing the spatial variability in peak season CO₂ exchange characteristics across the Arctic tundra using a light response curve parameterization.** *Biogeosciences* 11(17): 4897–4912.
- McVeigh, P., et al. (2014). **Meteorological and functional response partitioning to explain interannual variability of CO₂ exchange at an Irish Atlantic blanket bog.** *Agricultural and Forest Meteorology* 194: 8–19.
- Meijide, A., et al. (2014). **CO₂ and energy fluxes from an oil palm plantation in Sumatra, Indonesia.** EGU General Assembly.
- Meirelles, M., et al. (2015). **Carbon dioxide exchange in a tropical wet grassland.** *Wetlands Ecology and Management*. 23(5): 817–826.
- Mekhmandarov, Y., et al. (2015). **Application of the surface renewal technique in two types of screenhouses: Sensible heat flux estimates and turbulence characteristics.** *Agricultural and Forest Meteorology* 203: 229–242.
- Mengistu, M., et al. (2014). **The impact of taro (*Colocasia esculenta*) cultivation on the total evaporation of a *Cyperus latifolius* marsh.** *Hydrological Processes* 28(3): 620–627.
- Merbold, L., et al. (2014). **Greenhouse gas budget (CO₂, CH₄ and N₂O) of intensively managed grassland following restoration.** *Global Change Biology* 20(6): 1913–1928.
- Meyer, W. S., et al. (2015). **Evaporation of perennial semi-arid woodland in southeastern Australia is adapted for irregular but common dry periods.** *Hydrological Processes* 29(17): 3714–3726.
- Migliaccio, K. and W. Barclay Shoemaker (2014). **Estimation of urban subtropical bahiagrass (*Paspalum notatum*) evapotranspiration using crop coefficients and the eddy covariance method.** *Hydrological Processes* 28(15): 4487–4495.
- Miller, L. A., et al. (2015). **Methods for biogeochemical studies of sea ice: The state of the art, caveats, and recommendations.** *Elementa: Science of the Anthropocene* 3(1): 000038.
- Moat, B.I., et al. (2015). **Metadata for the WAGES instrumentation deployed on the RRS James Clark Ross between May 2010 and September 2013.** National Oceanography Centre.

- Monteiro, P., et al. (2014). **Estimation of energy balance components and evapotranspiration in soybean crop in southern Brazil using TM-Landsat 5 images.** *Bragantia* 73(1): 72–80.
- Monteiro, P., et al. **Evapotranspiration for soybean growing areas in central region of Rio Grande Do Sul State with data from TM sensor.** Conference paper, II Inovagri International Meeting.
- Moravek, A., et al. (2014). **Application of a GC-ECD for measurements of biosphere–atmosphere exchange fluxes of peroxyacetyl nitrate using the relaxed eddy accumulation and gradient method.** *Atmospheric Measurement Techniques* 7 (7): 2097–2119.
- Moravek, A., et al. (2015). **Influence of local air pollution on the deposition of peroxyacetyl nitrate to a nutrient-poor natural grassland ecosystem.** *Atmospheric Chemistry and Physics* 15 (2): 899–911.
- Morin, T., et al. (2014). **Environmental drivers of methane fluxes from an urban temperate wetland park.** *Journal of Geophysical Research: Biogeosciences* 119(11): 2188–2208.
- Morin, T., et al. (2014). **The seasonal and diurnal dynamics of methane flux at a created urban wetland.** *Ecological Engineering* 72: 74–83.
- Mouat, D., et al. (2014). **Carbon sequestration at United States Marine Corps installations west.** DTIC Document.
- Mutzner, R., et al. (2015). **Controls on the diurnal streamflow cycles in two subbasins of an alpine headwater catchment.** *Water Resources Research* 51(5): 3403–3418.
- Nakaji, T., et al. (2014). **Estimation of light-use efficiency through a combinational use of the photochemical reflectance index and vapor pressure deficit in an evergreen tropical rainforest at Pasoh, Peninsular Malaysia.** *Remote Sensing of Environment* 150: 82–92.
- Nakano, T. and M. Shinoda (2015). **Modeling gross primary production and ecosystem respiration in a semiarid grassland of Mongolia.** *Soil Science and Plant Nutrition* 61(1): 106–115.
- Ng, B. J., et al. (2014). **Carbon fluxes from an urban tropical grassland.** *Environmental Pollution* 203: 227–234.
- Novick, K. A., et al. (2015). **On the difference in the net ecosystem exchange of CO₂ between deciduous and evergreen forests in the southeastern United States.** *Global Change Biology*. 21: 827–842.
- Oechel, W. C., et al. (2014). **Annual patterns and budget of CO₂ flux in an Arctic tussock tundra ecosystem.** *Journal of Geophysical Research: Biogeosciences* 119(3): 323–339.
- Ohta, T., et al. (2014). **Effects of waterlogging on water and carbon dioxide fluxes and environmental variables in a Siberian larch forest, 1998–2011.** *Agricultural and Forest Meteorology* 188: 64–75.
- Oikawa, P. Y., et al. (2014). **Offsetting high water demands with high productivity: Sorghum as a biofuel crop in a high irradiance arid ecosystem.** GCB Bioenergy.
- Olchev, A., et al. (2015). **Response of CO₂ and H₂O fluxes of a mountainous tropical rain forest in equatorial Indonesia to El Niño events.** *Biogeosciences Discussions* 12: 4405–4431.
- Ono, K. and A. Maruyama (2015). **Development of an onsite computation scheme of eddy-covariance fluxes.** *Journal of Agricultural Meteorology*. Published online: 12 September 2015.
- Ouyang, Z., et al. (2014). **Disentangling the confounding effects of PAR and air temperature on net ecosystem exchange at multiple time scales.** *Ecological Complexity* 19: 46–58.
- Palatella, L., et al. (2014). **Towards a flux-partitioning procedure based on the direct use of high-frequency eddy-covariance data.** *Boundary-Layer Meteorology* 153(2): 327–337.
- Pardo, N., et al. (2014). **SEBS validation in a Spanish rotating crop.** *Agricultural and Forest Meteorology* 195: 132–142.
- Patil, M., et al. (2014). **Measurements of carbon dioxide and heat fluxes during monsoon-2011 season over rural site of India by eddy covariance technique.** *Journal of Earth System Science* 123(1): 177–185.
- Peichl, M., et al. (2014). **Carbon and greenhouse gas balances in an age sequence of temperate pine plantations.** *Biogeosciences* 11(19): 5399–5410.
- Pelletier, L., et al. (2014). **Carbon release from boreal peatland open water pools: Implication for the contemporary C exchange.** *Journal of Geophysical Research: Biogeosciences* 119(3): 207–222.
- Peng, Z. and J. Sun (2014). **Characteristics of the drag coefficient in the roughness sublayer over a complex urban surface.** *Boundary-Layer Meteorology* 153(3): 569–580.

- Personne, E., et al. (2015). **Investigating sources and sinks for ammonia exchanges between the atmosphere and a wheat canopy following slurry application with trailing hose.** *Agricultural and Forest Meteorology* 207: 11–23.
- Petrie, M. D. (2014). **Community ecology, climate change and ecohydrology in desert grassland and shrubland.** Dissertation: The University of New Mexico.
- Petrie, M., et al. (2014). **Grassland to shrubland state transitions enhance carbon sequestration in the northern Chihuahuan Desert.** *Global Change Biology* 21(3): 1226–1235.
- Petrie, M., et al. (2015). **The ecological role of small rainfall events in a desert grassland.** *Ecohydrology*. 1 January 2015. DOI: 10.1002/eco.1614.
- Petrone, R. M., et al. (2014). **Effects of harvesting and drought on CO₂ and H₂O fluxes in an aspen-dominated western boreal plain forest: early chronosequence recovery.** *Canadian Journal of Forest Research* 45(1): 87–100.
- Pinto, O. B., et al. (2014) **Energy balance and canopy conductance of a tropical semi-deciduous forest of the southern Amazon Basin.** *Water Resources Research* 44(3).
- Plake, D., et al. (2015). **Comparison of ozone deposition measured with the dynamic chamber and the eddy covariance method.** *Agricultural and Forest Meteorology* 206: 97–112.
- Podgrajsek, E., et al. (2014). **Comparison of floating chamber and eddy covariance measurements of lake greenhouse gas fluxes.** *Biogeosciences* 11 (15): 4225–4233.
- Podgrajsek, E., et al. (2015). **Diel cycle of lake-air CO₂ flux from a shallow lake and the impact of waterside convection on the transfer velocity.** *Journal of Geophysical Research: Biogeosciences* 120 (1): 29–38.
- Qi, Y., et al. (2015). **Using the Cross-Correlation Function to Evaluate the Quality of Eddy-Covariance Data.** *Boundary-layer Meteorology* 157 (2): 173–189.
- Quansah, E., et al. (2015). **Carbon dioxide fluxes from contrasting ecosystems in the Sudanian Savanna in West Africa.** *Carbon Balance and Management* 10(1): 1.
- Raddatz, R., et al. (2015). **A Simple Scheme for Estimating Turbulent Heat Flux over Landfast Arctic Sea Ice from Dry Snow to Advanced Melt.** *Boundary-Layer Meteorology* 155: 351–367.
- Rajan, N., et al. (2014). **Extreme drought effects on summer evapotranspiration and energy balance of a grassland in the Southern Great Plains.** *Ecohydrology*. First published online: 22 December 2014. DOI: 10.1002/eco.1574.
- Rajewski, D. A., et al. (2014). **Changes in fluxes of heat, H₂O, and CO₂ caused by a large wind farm.** *Agricultural and Forest Meteorology* 194: 175–187.
- Reed, D. E., et al. (2014). **Impact of mountain pine beetle induced mortality on forest carbon and water fluxes.** *Environmental Research Letters* 9(10): 105004.
- Ridler, M. E., et al. (2014). **Assimilation of SMOS - derived soil moisture in a fully integrated hydrological and soil-vegetation-atmosphere transfer model in Western Denmark.** *Water Resources Research* 50(11): 8962–8981.
- Riederer, M., et al. (2014). **Application of relaxed eddy accumulation (REA) on managed grassland.** *Atmospheric Measurement Techniques Discussions* 7: 4987–5026.
- Riederer, M., et al. (2014). **Net ecosystem CO₂ exchange measurements by the closed chamber method and the eddy covariance technique and their dependence on atmospheric conditions.** *Atmospheric Measurement Techniques* 7 (4): 1057–1064.
- Riederer, M., et al. (2014). **Prerequisites for application of hyperbolic relaxed eddy accumulation on managed grasslands and alternative net ecosystem exchange flux partitioning.** *Atmospheric Measurement Techniques* 7(12): 4237–4250.
- Riederer, M., et al. (2015). **Partitioning NEE for absolute C input into various ecosystem pools by combining results from eddy-covariance, atmospheric flux partitioning and ¹³C pulse labeling.** *Plant and Soil* 390(1–2): 61–76. ²
- Ringgaard, R., et al. (2014). **Partitioning forest evapotranspiration: Interception evaporation and the impact of canopy structure, local and regional advection.** *Journal of Hydrology* 517: 677–690.
- Rodrigues, T. R., et al. (2014). **Seasonal variation in energy balance and canopy conductance for a tropical savanna ecosystem of south central Mato Grosso, Brazil.** *Journal of Geophysical Research: Biogeosciences* 119(1): 1–13.
- Rutledge, S., et al. (2014). **CO₂ emissions following cultivation of a temperate permanent pasture.** *Agriculture, Ecosystems and Environment* 184: 21–33.

- Rutledge, S., et al. (2015). **Carbon balance of an intensively grazed temperate dairy pasture over four years.** *Agriculture, Ecosystems and Environment* 206: 10–20.
- Sahlée, E., et al. (2014). **Influence from surrounding land on the turbulence measurements above a lake.** *Boundary-Layer Meteorology* 150(2): 235–258.
- Sakowska, K., et al. (2014). **Monitoring of carbon dioxide fluxes in a subalpine grassland ecosystem of the Italian Alps using a multi-spectral sensor.** *Biogeosciences* 11(17): 4695–4712.
- Sakowska, K., et al. (2015). **WhiteRef: A new tower-based hyperspectral system for continuous reflectance measurements.** *Sensors* 15(1): 1088–1105.
- Sanchez-Mejia, Z. M. and S. A. Papuga (2014). **Observations of a two-layer soil moisture influence on surface energy dynamics and planetary boundary layer characteristics in a semiarid shrubland.** *Water Resources Research* 50(1): 306–317.
- Sandeep, A., et al. (2014). **Differences in atmospheric boundary-layer characteristics between wet and dry episodes of the Indian summer monsoon.** *Boundary-Layer Meteorology* 153(2): 217–236.
- Schäfer, K., et al. (2014). **Carbon dioxide fluxes of an urban tidal marsh in the Hudson-Raritan estuary.** *Journal of Geophysical Research: Biogeosciences* 119(11): 2065–2081.
- Scott, R. L., et al. (2014). **When vegetation change alters ecosystem water availability.** *Global Change Biology* 20(7): 2198–2210.
- Sen, L. and Z. Zhong (2014). **Improvement of surface flux calculation: A study based on measurements over alpine meadow in the eastern Tibet Plateau in summer.** *Chinese Physics B* 23(2): 029201.
- Senapati, N., et al. (2014). **Net carbon storage measured in a mowed and grazed temperate sown grassland shows potential for carbon sequestration under grazed system.** *Carbon Management* 5(2): 131–144.
- Serrano-Ortiz, P., et al. (2015). **Seasonality of net carbon exchanges of Mediterranean ecosystems across an altitudinal gradient.** *Journal of Arid Environments* 115: 1–9.
- Shafian, S. (2014). **Estimation of soil moisture status in the Texas High Plains using remote sensing.** Dissertation: Texas Tech University.
- Shang, L., et al. (2015). **Energy exchange of an alpine grassland on the eastern Qinghai-Tibetan Plateau.** *Science Bulletin*: 60(4) 435–446.
- Shimizu, T., et al. (2014). **Inter-annual variation of water vapor exchange measured by the bandpass eddy covariance method over a dry evergreen forest in the central Cambodia.** Proceedings of 10th International Workshop on Forest Watershed Environment Research in Cambodia.
- Shoemaker, W., et al. (2015). **Carbon exchange between the atmosphere and subtropical forested cypress and pine wetlands.** *Biogeosciences* 12(8): 2285–2300.
- Sievers, J., et al. (2015). **Estimating surface fluxes using eddy covariance and numerical optimization.** *Atmospheric Chemistry and Physics* 15(4): 2081–2103.
- Sievers, J., et al. (2015). **Winter observations of CO₂ exchange between sea-ice and the atmosphere in a coastal fjord environment.** *The Cryosphere Discussions* 9(1): 45–75.
- Skinner, R. H. and C. J. Dell (2015). **Comparing pasture C sequestration estimates from eddy covariance and soil cores.** *Agriculture, Ecosystems and Environment* 199: 52–57.
- Song, G., et al. (2014). **Micrometeorological flux measurements at a coastal site.** EGU General Assembly.
- Song, Q.H., et al. (2014). **Do the rubber plantations in tropical China act as large carbon sinks?** *Forest-Biogeosciences and Forestry* 7(1): 42–47.
- Song, W., et al. (2015). **Methane emissions from an alpine wetland on the Tibetan Plateau: Neglected but vital contribution of the nongrowing season.** *Journal of Geophysical Research: Biogeosciences* 120(8): 1475–1490.
- Sørensen, L. L., et al. (2014). **Parameterization of atmosphere–surface exchange of CO₂ over sea ice.** *Cryosphere* 8(3): 853–866.
- Souza, M. C., et al. (2014). **Estimation of gross primary production of the Amazon-Cerrado transitional forest by remote sensing techniques.** *Revista Brasileira de Meteorologia* 29(1): 1–12.
- Starkenbug, D., et al. (2015). **Temperature regimes and turbulent heat fluxes across a heterogeneous canopy in an Alaskan boreal forest.** *Journal of Geophysical Research: Atmospheres* 120(4): 1348–1360.

- Starr, G., et al. (2015). **Time series analysis of forest carbon dynamics: recovery of *Pinus palustris* physiology following a prescribed fire.** *New Forests* 46(1): 63–90.
- Sueki, S., et al. (2015). **Defoliation effects of *Diorhabda carinulata* on tamarisk evapotranspiration and groundwater levels.** *Ecohydrology*. 1 January 2015. DOI: 10.1002/eco.1604.
- Sugawara, H., et al. (2014). **How much does urban green cool town?** *Journal of Heat Island Institute International* 9: 2.
- Sun, J., et al. (2014). **Estimating Daytime Ecosystem Respiration to Improve Estimates of Gross Primary Production of a Temperate Forest.** *PLoS one* 9(11): e113512.
- Sun, K., et al. (2014). **On-road ammonia emissions characterized by mobile, open-path measurements.** *Environmental Science and Technology* 48(7): 3943–3950.
- Sun, K., et al. (2015). **Open-path eddy covariance measurements of ammonia fluxes from a beef cattle feedlot.** *Agricultural and Forest Meteorology*, 213: 193–202.
- Sun, K., et al. (2015). **Quantifying the Influence of Random Errors in Turbulence Measurements on Scalar Similarity in the Atmospheric Surface Layer.** *Boundary-layer Meteorology* 157(1): 61–80.
- Sun, K., et al. (2015). **Validation of TES ammonia observations at the single pixel scale in the San Joaquin Valley during DISCOVER-AQ.** *Journal of Geophysical Research: Atmospheres* 120(10): 5140–5154.
- Sun, W., et al. (2014). **Ablation modeling and surface energy budget in the ablation zone of Laohugou glacier No. 12, western Qilian mountains, China.** *Annals of Glaciology* 55(66): 111–120.
- Tagesson, T., et al. (2015). **Dynamics in carbon exchange fluxes for a grazed semi-arid savanna ecosystem in West Africa.** *Agriculture, Ecosystems and Environment* 205: 15–24.
- Tan, Z. H., et al. (2015). **Interannual and seasonal variability of water use efficiency in a tropical rainforest: Results from a 9 year eddy flux time series.** *Journal of Geophysical Research: Atmospheres* 120(2): 464–479.
- Tang, R., et al. (2015). **On the feasibility of temporally upscaling instantaneous evapotranspiration using weather forecast information.** *International Journal of Remote Sensing*: First published online: 7 April 2015. DOI:10.1080/01431161.2015.1029597.
- Tang, Y., et al. (2014). **Interannual variation of the Bowen ratio in a subtropical coniferous plantation in Southeast China, 2003–2012.** *PLoS one* 9(2): e88267.
- Tang, Y., et al. (2014). **The limiting effect of deep soilwater on evapotranspiration of a subtropical coniferous plantation subjected to seasonal drought.** *Advances in Atmospheric Sciences* 31(2): 385–395.
- Tao, L., et al. (2015). **Low-power, open-path mobile sensing platform for high-resolution measurements of greenhouse gases and air pollutants.** *Applied Physics B* 119(1): 153–164.
- Taylor, N., et al. (2014). **Crop coefficient approaches based on fixed estimates of leaf resistance are not appropriate for estimating water use of citrus.** *Irrigation Science* 33: 153–166.
- Thorp, R. (2014). **Observations of heat, water vapor and carbon dioxide exchanges over a living roof using eddy covariance.** Thesis: San Francisco State University.
- Tian, S., et al. (2015). **Testing DRAINMOD-Forest for predicting evapotranspiration in a mid-rotation pine plantation.** *Forest Ecology and Management*.
- Timm, A. U., et al. (2014). **Energy Partitioning and Evapotranspiration over a Rice Paddy in Southern Brazil.** *Journal of Hydrometeorology* 15(5): 1975–1988.
- Tokoro, T., et al. (2014). **Net uptake of atmospheric CO₂ by coastal submerged aquatic vegetation.** *Global Change Biology* 20(6):1873–1884.
- Tong, X., et al. (2014). **Biophysical Controls on Light Response of Net CO₂ Exchange in a Winter Wheat Field in the North China Plain.** *PLoS one* 9 (2): e89469.
- Tong, X., et al. (2014). **Ecosystem water use efficiency in a warm-temperate mixed plantation in the North China.** *Journal of Hydrology* 512: 221–228.
- Torn, M. and D. B. N. Raz-Yaseef (2014). **Eddy-Covariance and auxiliary measurements, NGEE-Barrow, 2012–2013.** Next Generation Ecosystems Experiment-Arctic, Oak Ridge National Laboratory (ORNL), Oak Ridge, TN (US).
- Trant, J. S. (2014). **Effects of thinning on carbon dynamics in a temperate coniferous forest.** Thesis: McMaster University.
- Ueyama, M., et al. (2014). **Autumn warming reduces the CO₂ sink of a black spruce forest in interior Alaska based on a nine-year eddy covariance measurement.** *Global Change Biology* 20(4):1161–1173.

- Ueyama, M., et al. (2014). **Inferring methane fluxes at a larch forest using Lagrangian, Eulerian, and hybrid inverse models.** *Journal of Geophysical Research: Biogeosciences* 119(10): 2018–2031.
- ÜNLÜ, M., et al. (2014). **Irrigation scheduling of grapefruit trees in a Mediterranean environment throughout evaluation of plant water status and evapotranspiration.** *Turkish Journal of Agriculture and Forestry* 38(6): 908–915.
- van Asperen, H., et al. (2015). **The role of photo-and thermal degradation for CO₂ and CO fluxes in an arid ecosystem.** *Biogeosciences Discussions* 12(3): 2429–2457.
- van de Boer, A., et al. (2014). **Estimation of the refractive index structure parameter from single-level daytime routine weather data.** *Applied Optics* 53: 5944–5960.
- Vásquez, V., et al. (2015). **Integrating lysimeter drainage and eddy covariance flux measurements in a groundwater recharge model.** *Hydrological Sciences Journal*. 60(9): 1520–1537.
- Velasco, E., et al. (2014). **Sources and sinks of carbon dioxide in a neighborhood of Mexico City.** *Atmospheric Environment* 97: 226–238.
- Velluet, C., et al. (2014). **Building a field-and model-based climatology of local water and energy cycles in the cultivated Sahel—Annual budgets and seasonality.** *Hydrology and Earth System Sciences* 18(12): 5001–5024.
- Verkerke, J. L., et al. (2014). **Remote sensing of CO₂ leakage from geologic sequestration projects.** *International Journal of Applied Earth Observation and Geoinformation* 31: 67–77.
- Vourlitis, G. L., et al. (2014). **Variations in evapotranspiration and climate for an Amazonian semi-deciduous forest over seasonal, annual, and El Niño cycles.** *International Journal of Biometeorology* 59(2): 217–230.
- Wagle, P. and V. G. Kakani (2014). **Confounding effects of soil moisture on the relationship between ecosystem respiration and soil temperature in switchgrass.** *BioEnergy Research* 7(3): 789–798.
- Wagle, P. and V. G. Kakani (2014). **Environmental control of daytime net ecosystem exchange of carbon dioxide in switchgrass.** *Agriculture, Ecosystems and Environment* 186: 170–177.
- Wagle, P. and V. G. Kakani (2014). **Growing season variability in evapotranspiration, ecosystem water use efficiency, and energy partitioning in switchgrass.** *Ecohydrology* 7(1): 64–72.
- Wagle, P. and V. G. Kakani (2014). **Seasonal variability in net ecosystem carbon dioxide exchange over a young Switchgrass stand.** *GCB Bioenergy* 6(4): 339–350.
- Wagle, P., et al. (2015). **Net ecosystem carbon dioxide exchange of dedicated bioenergy feedstocks: Switchgrass and high biomass sorghum.** *Agricultural and Forest Meteorology* 207: 107–116.
- Wang, J., et al. (2014). **Assessment of Uncertainties in Eddy Covariance Flux Measurement Based on Intensive Flux Matrix of HiWATER-MUSOEXE.** *Geoscience and Remote Sensing Letters, IEEE* 12(2): 259–263
- Wang, L., et al. (2015). **Contrasting characteristics of the surface energy balance between the urban and rural areas of Beijing.** *Advances in Atmospheric Sciences* 32(4): 505–514.
- Wang, L., et al. (2014). **Turbulent transport of momentum and scalars above an urban canopy.** *Boundary-Layer Meteorology* 150(3): 485–511.
- Wang, L., et al. (June 10, 2015). **Water and carbon dioxide fluxes over an alpine meadow in southwest China and the impact of a spring drought event.** *International Journal of Biometeorology*. First published online: 10 June 2015.
- Wang, X., et al. (2014). **Wind regimes above and below a temperate deciduous forest canopy in complex terrain: Interactions between slope and valley winds.** *Atmosphere* 6(1): 60–87.
- Wang, Y., et al. (2014). **Observed and simulated features of the CO₂ diurnal cycle in the boundary layer at Beijing and Hefei, China.** *Chinese Science Bulletin* 59(14): 1529–1535.
- Wang, Y., et al. (2015). **Structures of convection and turbulent kinetic energy in boundary layer over the southeastern edge of the Tibetan Plateau.** *Science China Earth Sciences* 58(7): 1198–1209.
- Wang, Z., et al. (2015). **Quantifying changes in multiple ecosystem services during 1992–2012 in the Sanjiang Plain of China.** *Science of The Total Environment* 514: 119–130.
- Ward, H., et al. (2014). **Multi-Scale Sensible Heat Fluxes in the Suburban Environment from Large-Aperture Scintillometry and Eddy Covariance.** *Boundary-Layer Meteorology* 152(1): 65–89.
- Ward, H., et al. (2015). **Effects of urban density on carbon dioxide exchanges: Observations of dense urban, suburban and woodland areas of southern England.** *Environmental Pollution* 198: 186–200.

- Ward, H., et al. (2015). **Infrared and millimetre-wave scintillometry in the suburban environment–Part 2: Large-area sensible and latent heat fluxes.** *Atmospheric Measurement Techniques* 8(3): 1407–1424.
- Ward, H., et al. (2015). **Infrared and millimetre-wave scintillometry in the suburban environment–Part 1: Structure parameters.** *Atmospheric Measurement Techniques* 8(3): 1385–1405.
- Wei, Z., et al. (2015). **Partitioning of evapotranspiration using high frequency water vapor isotopic measurement over a rice paddy field.** *Water Resources Research*. 51(5): 3716–372.
- Wharton, S., et al. (2014). **Role of surface energy exchange for simulating wind turbine Inflow: A case study in the southern great plains, USA.** *Atmosphere* 6(1): 21–49.
- Wickström, S. (2015). **Turbulent CO₂ exchange and surface energy balance over melting fjord-ice in North-East Greenland.** Masters thesis: University of Helsinki Department of Physics.
- Wildmann, N., et al. (2014). **An inverse-modelling approach for frequency response correction of capacitive humidity sensors in ABL research with small remotely piloted aircraft (RPA).** *Atmospheric Measurement Techniques* 7(9): 3059–3069.
- Wilson, T., et al. (2014). **Leaf litter water content and soil surface CO₂ fluxes in a deciduous forest.** *Agricultural and Forest Meteorology* 192: 42–50.
- Xiao, W., et al. (2014). **A flux-gradient system for simultaneous measurement of the CH₄, CO₂, and H₂O fluxes at a lake–air interface.** *Environmental Science and Technology* 48(24): 14490–14498.
- Xiao-Feng, G. (2014). **Effects of coordinate rotation on turbulent flux measurements during wintertime haze pollution in Beijing, China.** *Atmospheric and Oceanic Science Letters*: 8(2): 67–71.
- Xie, J., et al. (2014). **Long-term variability in the water budget and its controls in an oak dominated temperate forest.** *Hydrological Processes* 28(25): 6054–6066.
- Xie, X., et al. (2014). **Dependence of coastal wetland ecosystem respiration on temperature and tides: a temporal perspective.** *Biogeosciences* 11 (3): 539–545.
- Xu, L., et al. (2014). **Impact of changes in barometric pressure on landfill methane emission.** *Global Biogeochemical Cycles* 28(7): 679–695.
- Xu, M., et al. (2014). **Effects of climatic factors and ecosystem responses on the inter-annual variability of evapotranspiration in a coniferous plantation in subtropical China.** *PloS one* 9(1): e85593.
- Xue, B. L., et al. (2014). **Canopy conductance in a two-storey Siberian boreal larch forest, Russia.** *Hydrological Processes* 29(6): 1017–1026.
- Yan, G., et al. (2014). **Seasonal changes of energy fluxes of a subtropical mangrove forest in Zhangjiang Estuary, China.** *Journal of Chemical and Pharmaceutical Research* 6(1): 363–368.
- Yang, F., et al. (2014). **Evapotranspiration Measurement and Crop Coefficient Estimation over a Spring Wheat Farmland Ecosystem in the Loess Plateau.** *PloS one* 9(6): e100031.
- Yao, J., et al. (2014). **The surface energy budget on the debris-covered Koxkar Glacier in China.** *Environmental Earth Sciences* 72(11): 4503–4510.
- Ye, X., et al. (2015). **The turbulent structure and transport in fog layers observed over the Tianjin area.** *Atmospheric Research* 153: 217–234.
- Yu, Q., et al. (2014). **Narrowband bio-indicator monitoring of temperate forest carbon fluxes in northeastern China.** *Remote Sensing* 6(9): 8986–9013.
- Yuan, W., et al. (2015). **Uncertainty in simulating gross primary production of cropland ecosystem from satellite-based models.** *Agricultural and Forest Meteorology* 207: 48–57.
- Zhang, G., et al. (2014). **Analysis of the variability of canopy resistance over a desert steppe site in Inner Mongolia, China.** *Advances in Atmospheric Sciences* 31(3): 681–692.
- Zhang, G., et al. (2014). **Multi-scale decomposition of turbulent fluxes above a forest canopy.** *Agricultural and forest meteorology* 186: 48–63.
- Zhang, H. and H. Zhang (2015). **Comparison of turbulent sensible heat flux determined by large-aperture scintillometer and eddy covariance over urban and suburban areas.** *Boundary-Layer Meteorology* 154(1): 119–136.
- Zhang, L. (2014). **Quantification of fugitive emissions from a biosolids lagoon.** Master thesis: University of Alberta.
- Zhang, L., et al. (2015). **Diurnal and seasonal variations in carbon dioxide exchange in ecosystems in the zhangye oasis area, Northwest China.** *PloS one* 10(3): e0120660.

- Zhang, Q., et al. (2015). **Characteristics of Dew Formation and Distribution, and Its Contribution to the Surface Water Budget in a Semi-arid Region in China.** *Boundary-layer Meteorology* 154 (2): 317–331.
- Zhang, Q. and H. Liu (2014). **Seasonal changes in physical processes controlling evaporation over inland water.** *Journal of Geophysical Research: Atmospheres* 119(16): 9779–9792.
- Zhang, T., et al. (2014). **Lagged climatic effects on carbon fluxes over three grassland ecosystems in China.** *Journal of Plant Ecology* 8(3).
- Zhao, J., et al. (2014). **Spatial representativeness of eddy covariance measurements using footprint analysis in arid areas.** *Environmental earth sciences* 71(4): 1691–1697.
- Zhao, P., et al. (2015). **Comparison of dual crop coefficient method and Shuttleworth-Wallace model in evapotranspiration partitioning in a vineyard of northwest China.** *Agricultural Water Management* 160: 41–56.
- Zhao, P. (2014). **Ecosystem-atmosphere exchange of carbon dioxide and water vapour in typical East-Asian croplands.** Thesis, University of Bayreuth.
- Zheng, D., et al. (2014). **Assessment of roughness length schemes implemented within the Noah land surface model for high-altitude regions.** *Journal of Hydrometeorology* 15(3): 921–937.
- Zheng, H., et al. (2014). **Hysteresis Responses of Evapotranspiration to Meteorological Factors at a Diel Timescale: Patterns and Causes.** *PLoS one* 9(6): e98857.
- Zhou, J., et al. (2014). **Water-use efficiency of a poplar plantation in Northern China.** *Journal of Forest Research* 19(6): 483–492.
- Zhou, X., et al. (2014). **Comparison of ET estimations by the three-temperature model, SEBAL model and eddy covariance observations.** *Journal of Hydrology* 519: 769–776.
- Zhu, G., et al. (2014). **Modelling evapotranspiration in an alpine grassland ecosystem on Qinghai-Tibetan plateau.** *Hydrological Processes* 28(3): 610–619.
- Zhu, G., et al. (2014). **Simultaneously assimilating multivariate data sets into the two-source evapotranspiration model by Bayesian approach: application to spring maize in an arid region of northwestern China.** *Geoscientific Model Development* 7(4): 1467–1482.
- Zhu, W., et al. (2014). **Air-surface exchange of Hg⁰ measured by collocated micrometeorological and enclosure methods—Part 1: Data comparability and method characteristics.** *Atmospheric Chemistry and Physics Discussions* 14 (16): 22273–22319.
- Zhu, X., et al. (2014). **Seasonal dynamics of water use efficiency of typical forest and grassland ecosystems in China.** *Journal of Forest Research* 19 (1): 70–76.
- Zhu, X.-J., et al. (2015). **Spatiotemporal variations of T/ET (the ratio of transpiration to evapotranspiration) in three forests of Eastern China.** *Ecological Indicators* 52: 411–421.
- Zhu, Z., et al. (2014). **Diurnal variation of ozone flux over corn field in Northwestern Shandong Plain of China.** *Science China Earth Sciences* 57(3): 503–511.
- Zielis, S., et al. (2014). **NEP of a Swiss subalpine forest is significantly driven not only by current but also by previous year's weather.** *Biogeosciences* 11(6): 1627–1635.
- Zona, D., et al. (2014). **Delayed responses of an Arctic ecosystem to an extreme summer: impacts on net ecosystem exchange and vegetation functioning.** *Biogeosciences* 11(20): 5877–5888.
- 강남구 and 윤주열 (2015). "개회로 파장 변조 분광법과 에디 공분산 방법으로 논에서 관측된 CH₄ 플럭스 자료의 보정." *한국농림기상학회지* 17(1): 15–24.
- 심교문, et al. "김제벼-보리 이모작 논에서 벼 재배기간 동안의 순생태계 CO₂ 교환량에 대한 환경요인 분석."
- 刘斌, et al. (2015). "水层厚度变化对稻田能量分配的影响." *生态学杂志* 34(3): 695–702.
- 刘玉莉, et al. (2014). "安吉毛竹林水汽通量变化特征及其与环境因子的关系." *生态学报* 34(17): 4900–4909.
- 刘郁珏, et al. (2014). "北京城市通量足迹及源区分布特征分析." *大气科学* 38(6): 1044–1054.
- 孙成, et al. (2014). "我国亚热带毛竹林生长季能量通量过程及闭合度分析." *热带亚热带植物学报* 22 (1): 38–44.
- 张摇蕾 (2014). "张掖灌区玉米农田生态系统 CO₂ 通量的变化规律和环境响应." *Ecology* 33(7): 1722–1728.
- 李炬 and 窦军霞 (2014). "北京城市气象观测试验进展." *气象科技进展* 1: 009.
- 李祎君 (2015). "雨养玉米农田水热交换的环境控制机理研究." *气候与环境研究* 20(1): 71–79.

- 王江涛, et al. (2014). "崇明东滩滨海围垦湿地 CO₂ 通量贡献区分析." *生态与农村环境学报* 30(5): 588–594.
- 王记明, et al. (2014). "青海湖高寒湿地生态系统夏季 CO₂ 通量日变化及其影响因子研究." *生态与农村环境学报* 30(3): 317–323.
- 贺俊杰 (2014). "典型草原 CO₂ 通量变化特征对环境因子的响应." *中国农学通报* 30(34): 107–111.
- 贺俊杰 (2014). "典型草原区草地潜热通量特征." *中国农学通报* 30(26): 195–199.
- 贾志军, et al. (2014). "三江平原大豆田蒸散量模拟." *生态与农村环境学报* 30(1): 50–56.
- 邢庆会, et al. (2014). "黄河口潮间盐沼湿地生长季净生态系统 CO₂ 交换特征及其影响因素."
- 邹陈, et al. (2014). "乌兰乌苏绿洲棉田生态系统水汽和二氧化碳通量研究." *中国农学通报* 30(15): 120–125.
- 郭阳, et al. (2015). "均匀裸土地表土壤热通量计算方法对比及对能量闭合的影响." *气候与环境研究* 20(2): 177–187.
- 马金龙, et al. (2015). "干旱区绿洲膜下滴灌棉田蒸散过程." *生态学杂志* 34(4): 974–981.
- Anderson, R. G., et al. (2015). **Long-rotation sugarcane in Hawaii sustains high carbon accumulation and radiation use efficiency in 2nd year of growth.** *Agriculture, Ecosystems and Environment* 199: 216–224.
- Artigas, F., et al. (2015). **Long term carbon storage potential and CO₂ sink strength of a restored salt marsh in New Jersey.** *Agricultural and Forest Meteorology* 200: 313–321.
- Begashaw, I. (2014). **A New Tool for Automated Data Collection and Complete On-Site Flux Data Processing for Eddy Covariance Measurements.** AGU Fall Meeting.
- Brandt, P., et al. (2015). **On the role of circulation and mixing in the ventilation of oxygen minimum zones with a focus on the eastern tropical North Atlantic.** *Biogeosciences* 12: 489–512.
- Buckley, S. M., et al. (2014). **Variations in carbon dioxide fluxes within a city landscape: Identifying a vehicular influence.** *Urban Ecosystems*. First published online: 1 February 2014.
- Campbell, D. I., et al., (2015). **Variations in CO₂ exchange for dairy farms with year-round rotational grazing on drained peatlands.** *Agriculture, Ecosystems and Environment* 202: 68–78.
- Chamberlain, S. D., et al. (2015). **Underlying Ecosystem Emissions Exceed Cattle-Emitted Methane from Subtropical Lowland Pastures.** *Ecosystems* 18(6): 933–945.
- Coates, T., et al. (2014). **A year In the life: The challenges of long term methane flux measurements on a cattle-grazed landscape.** *Optical Instrumentation for Energy and Environmental Applications*, Optical Society of America.
- Dangwala, N., et al. (2015). **Monitoring of water stress in wheat using multispectral indices derived from Landsat-TM.** *Geocarto International*. Published online: 1 September 2015.
- El-Madany, T.-S., et al. (2014). **Low-level jets and above canopy drainage as causes for turbulent exchange in the nocturnal boundary layer.** *Biogeosciences Discussions* 11(3): 4695–4727.
- Emmerton, C. A., et al. (2015). **Net ecosystem exchange of CO₂ with rapidly changing high Arctic landscapes.** *Global Change Biology*. First published online: 17 August 2015. DOI: 10.1111/gcb.13064.

EddyPro® Eddy Covariance Processing Software

- Ahongshangbam, J. (2014). **Carbon flux monitoring and modelling in terai central forest division, Haldwani.** Thesis: Andhra University, Visakhapatnam.
- Alberto, M. C. R., et al. (2014). **Actual evapotranspiration and dual crop coefficients for dry-seeded rice and hybrid maize grown with overhead sprinkler irrigation.** *Agricultural Water Management* 136(4): 1–12.
- Alberto, M. C. R., et al. (2014). **Measuring methane flux from irrigated rice fields by eddy covariance method using open-path gas analyzer.** *Field Crops Research* 160: 12–21.
- Anderson, R. G., et al. (2015). **Divergence of actual and reference evapotranspiration observations for irrigated sugarcane with windy tropical conditions.** *Hydrology and Earth System Sciences* 19(1): 583–600.

- Emmerton, C. A. (2015). **The net exchange of carbon greenhouse gases with high Arctic terrestrial and aquatic ecosystems.** Doctoral Thesis: Department of Biological Sciences, University of Alberta.
- Falocchi, M., et al. (2015). **Turbulence measurements in an Alpine valley: The CividatEX Experiment case.** EGU General Assembly.
- Fleischer, E., et al. (2015). **Summer evapotranspiration in western Siberia: a comparison between eddy covariance and Penman method formulations.** *Hydrological Processes* 29(20): 4498–4513.
- Fratini, G., D. K. McDermitt, and D. Papale (2014). **Eddy-covariance flux errors due to biases in gas concentration measurements: Origins, quantification and correction.** *Biogeosciences* 11 (4): 1037–1051.
- Fratini, G., and Mauder, M. (2014). **Towards a consistent eddy-covariance processing: an intercomparison of EddyPro and TK3.** *Atmospheric Measurement Techniques Discussions* 7 (3): 2107–2126.
- Gao, G., et al. (2015). **Comparison of three evapotranspiration models with eddy covariance measurements for a *Populus euphratica* Oliv. forest in an arid region of northwestern China.** *Journal of Arid Land*. First published online: 30 August 2015.
- Gong, T. T., et al. (2015). **Characteristics of surface energy fluxes over a sparse shrubland ecosystem in the farming-pastoral zone of the Loess Plateau, Northwest China.** *Chinese Journal of Applied Ecology* 26(6): 1625–1633.
- Gong, T. T., et al. (2014). **Effects of vegetation change on evapotranspiration in a semiarid shrubland of the Loess Plateau, China.** *Hydrology and Earth System Sciences Discussions* 11(12): 13571–13605.
- Goodrich, J. P., et al. (2015). **High vapor pressure deficit constrains GPP and the light response of NEE at a Southern Hemisphere bog.** *Agricultural and Forest Meteorology* 203: 54–63.
- Goodrich, J. P., et al. (2015). **Overriding control of methane flux temporal variability by water table dynamics in a Southern Hemisphere, raised bog.** *Journal of Geophysical Research: Biogeosciences* 120(5): 819–831.
- Herbst, M., et al. (2015). **Differences in carbon uptake and water use between a managed and an unmanaged beech forest in central Germany.** *Forest Ecology and Management* 355: 101–108.
- Hoffmann, H., et al. (2015). **Estimating evapotranspiration with thermal UAV data and two source energy balance models.** *Hydrology and Earth System Sciences Discussions* 12: 7469–7502.
- Hommeltenberg, J., et al. (2014). **Ecosystem scale methane fluxes in a natural temperate bog-pine forest in southern Germany.** *Agricultural and Forest Meteorology* 198: 273–284.
- Huang, H., et al. (2015). **Atmospheric boundary layer structure and turbulence during sea fog on the southern China coast.** *Monthly Weather Review* 143: 1907–1923
- Huang, H., et al. (2014). **Nitrous oxide emissions from a commercial cornfield (*Zea mays*) measured using the eddy-covariance technique.** *Atmospheric Chemistry and Physics Discussions* 14 (14): 20417–20460.
- Huo, Q., et al. (2015). **Effects of surface source/sink distributions on the flux-gradient similarity in the unstable surface layer.** *Theoretical and Applied Climatology* 119: 313–322.
- Hurkuck, M., et al. (2014). **Determination of atmospheric nitrogen deposition to a semi-natural peat bog site in an intensively managed agricultural landscape.** *Atmospheric Environment* 97: 296–309.
- Ikawa, H., et al. (2015). **Understory CO₂, sensible heat, and latent heat fluxes in a black spruce forest in interior Alaska.** *Agricultural and Forest Meteorology* 214-215: 80-90.
- Jha, C. S., et al. (2014). **Eddy covariance based methane flux in Sundarbans mangroves, India.** *Journal of Earth System Science : Published by the Indian Academy of Sciences* 123(5): 1089–1096.
- Ingwersen, J., et al. (2015). **On the use of the post-closure methods uncertainty band to evaluate the performance of land surface models against eddy covariance flux data.** *Biogeosciences* 12: 2311–2326.
- Isabelle, P., et al., (2015). **Applicability of the Bulk-Transfer Approach to Estimate Evapotranspiration from Boreal Peatlands.** *Journal of Hydrometeorology* 16: 1521–1539.
- Jia, X., et al. (2014). **Biophysical controls on net ecosystem CO₂ exchange over a semiarid shrubland in northwest China.** *Biogeosciences* 11 (17): 4679–4693.

- Jiemin, W., et al. (2015). **Assessment of Uncertainties in Eddy Covariance Flux Measurement Based on Intensive Flux Matrix of HiWATER-MUSOEXE.** *IEEE Geoscience and Remote Sensing Letters* 12(2): 259–263.
- Kiese, R., et al., (2015). **Biosphere-Atmosphere Exchange of CO₂ and CH₄ over Natural and Drained Bog Forest Ecosystems in Southern Germany.** Dissertation: Universitätsbibliothek der TU München.
- Lewicki, J. L. and G. E. Hilley, (2014). **Multi-scale observations of the variability of magmatic CO₂ emissions, Mammoth Mountain, CA, USA.** *Journal of Volcanology and Geothermal Research*, 284: 1–15.
- Li, Z., et al., (2015). **Long-term energy flux and radiation balance observations over Lake Ngoring, Tibetan Plateau.** *Atmospheric Research* 155: 13–25.
- Li, Z., et al. (2015). **Turbulent transfer coefficient and roughness length in a high-altitude lake, Tibetan Plateau.** *Theoretical and Applied Climatology*. 17 April 2015.
- Liu, H. Z., et al. (2015). **Eddy covariance measurements of water vapor and CO₂ fluxes above the Erhai Lake.** *Science China Earth Sciences* 58(3): 317–328.
- Ma, N., et al. (2015). **Modeling actual evapotranspiration with routine meteorological variables in the data-scarce region of the Tibetan Plateau: Comparisons and implications.** *Journal of Geophysical Research: Biogeosciences* 120(8): 1638–1657.
- Macáľková, L. et al. (2015). **Flux footprints in different ecosystems.** Global Change: A Complex Challenge, Conference Proceedings. 54–60.
- Masseroni, D., et al., (2014). **Surface energy flux measurements in a flooded and an aerobic rice field using a single eddy-covariance system.** *Paddy and Water Environment*: 13(4): 405–424.
- Mazirov, I., et al. (2015). **Agricultural crops and soil treatment impacts on the daily and seasonal dynamics of CO₂ fluxes in the field agro-ecosystems at the Central region of Russia.** EGU General Assembly.
- Meshalkina, J., et al. (2015). **Central Russia agroecosystem monitoring with CO₂ fluxes analysis by eddy covariance method.** *Eurasian Journal of Soil Science* 4(3): 211.
- Mitchell, S. R., et al. (2015) **Land-atmosphere carbon and water flux relationships to vapor pressure deficit, soil moisture, and stream flow.** *Agricultural and Forest Meteorology* 208: 108–117.
- Nogu, Liisi. (2014). **The effects of site preparation on carbon fluxes at two clear-cuts in southern Sweden.** Master degree thesis: Department of Physical Geography and Ecosystems Science, Lund University.
- Owen, N. A., et al. (2015). **Eddy covariance captures four-phase crassulacean acid metabolism (CAM) gas exchange signature in Agave.** *Plant, Cell and Environment*. First published online: 8 September 2015. DOI: 10.1111/pce.12610.
- Rajan, N., et al. (2014). **Extreme drought effects on summer evapotranspiration and energy balance of a grassland in the Southern Great Plains.** *Ecohydrology*. First published online: 22 December 2014. DOI: 10.1002/eco.1574.
- Roll, A., et al. (2015). **Transpiration in an oil palm landscape: Effects of palm age.** *Biogeosciences Discussions*: 12: 9209–9242.
- Sakagami, Y., et al. (2015). **Effects of turbulence, wind shear, wind veer, and atmospheric stability on power performance: a case study in Brazil.** Santa Catarina Federal Institute. Florianópolis, Brazil.
- Shang, L., et al. (2015). **Energy exchange of an alpine grassland on the eastern Qinghai-Tibetan Plateau.** *Science Bulletin* 60(4): 435–446.
- Shoemaker, W. B., et al. (2015). **Carbon exchange between the atmosphere and subtropical forested cypress and pine wetlands.** *Biogeosciences* 12(8): 2285–2300.
- Song, W., et al. (2015). **Methane emissions from an alpine wetland on the Tibetan Plateau: Neglected but vital contribution of the nongrowing season.** *Journal of Geophysical Research: Biogeosciences* 120(8): 1475–1490.
- Sueki, S., et al. (2015). **Defoliation effects of *Diorhabda carinulata* on tamarisk evapotranspiration and groundwater levels.** *Ecohydrology*. First published online: 6 February 2015. DOI: 10.1002/eco.1604.
- Sun, K., et al. (2015). **Open-path eddy covariance measurements of ammonia fluxes from a beef cattle feedlot.** *Agricultural and Forest Meteorology* 213: 193–202.
- Tagesson, T., et al. (2015). **Deriving seasonal dynamics in ecosystem properties of semi-arid savanna grasslands from in situ-based hyperspectral reflectance.** *Biogeosciences* 12: 4621–4635.

- Tagesson, T., et al. (2015). **Dynamics in carbon exchange fluxes for a grazed semi-arid savanna ecosystem in West Africa.** *Agriculture, Ecosystems and Environment* 205: 15–24.
- Wagle, P. and V. G. Kakani (2014). **Environmental control of daytime net ecosystem exchange of carbon dioxide in switchgrass.** *Agriculture, Ecosystems and Environment* 186: 170–177.
- Wagle, P., et al. (2015). **Net Ecosystem Carbon Dioxide Exchange of Dedicated Bioenergy Feedstocks: Switchgrass and High Biomass Sorghum.** *Agricultural and Forest Meteorology* 207: 107–116.
- Wang, L., et al. (2015). **Water and carbon dioxide fluxes over an alpine meadow in southwest China and the impact of a spring drought event.** *International Journal of Biometeorology*. First published online: 10 June 2015.
- Wang, Y., et al. (2015). **Structures of convection and turbulent kinetic energy in boundary layer over the southeastern edge of the Tibetan Plateau.** *Science China Earth Sciences* 58(7): 1198–1209.
- Ward, H. C., et al. (2015). **Effects of urban density on carbon dioxide exchanges: Observations of dense urban, suburban and woodland areas of southern England.** *Environmental Pollution*: 198: 186–200.
- Ward, H.C., et al. (2015). **Infrared and millimetre-wave scintillometry in the suburban environment - Part 2: Large-area sensible and latent heat fluxes.** *Atmospheric Measurement Techniques* 8: 1407–1424.
- Yaroslavtsev, A., et al. (2015). **Climate determined differences in carbon dioxide fluxes dynamics between two comparable agroecosystems of Central Russia.** EGU General Assembly.
- Yee, M. S., et al. (2015). **A comparison of optical and microwave scintillometers with eddy covariance derived surface heat fluxes.** *Agricultural and Forest Meteorology* 213(5): 226–239.
- Zhang, He., and Hongsheng Zhang (2015). **Comparison of Turbulent Sensible Heat Flux Determined by Large-Aperture Scintillometer and Eddy Covariance over Urban and Suburban Areas.** *Boundary-layer Meteorology* 154(1): 119–136.
- Zhang, L. (2014). **Quantification of fugitive emissions from a biosolids lagoon.** Thesis: University of Alberta.
- Zhu, W., et al. (2015). **Mercury vapor air–surface exchange measured by collocated micro-meteorological and enclosure methods – Part II: Bias and uncertainty analysis.** *Atmospheric Chemistry and Physics* 15(10): 5359–5376.
- Zhu, Z., et al. (2015). **The effects of different calibration and frequency response correction methods on eddy covariance ozone flux measured with a dry chemiluminescence analyzer.** *Agricultural and Forest Meteorology* 213: 114–125.
- Zona, D., et al. (2014). **Environmental controls on ozone fluxes in a poplar plantation in Western Europe.** *Environmental Pollution* 184(5): 201–210.

Please contact us with any corrections or potential additions to this list.
Tell us about your research by visiting www.licor.com/case-study.

LI-COR Biosciences

Global Headquarters

4647 Superior Street • Lincoln, Nebraska 68504 • Phone: +1-402-467-3576 • Toll free: 800-447-3576
Fax: +1-402-467-2819 envsales@licor.com • envsupport@licor.com • www.licor.com/env

The LI-COR logo consists of the words "LI-COR" in a bold, italicized, sans-serif font. The "LI" is in a larger font size than "COR". A registered trademark symbol (®) is located at the top right of the "OR".

Regional Offices

LI-COR GmbH, Germany

Serving Andorra, Albania, Cyprus, Estonia, Germany, Iceland, Latvia, Lithuania, Liechtenstein, Malta, Moldova, Monaco, San Marino, Ukraine, and Vatican City.
LI-COR Biosciences GmbH • Siemensstraße 25A • 61352 Bad Homburg • Germany
Phone: +49 (0) 6172 17 17 771 • Fax: +49 (0) 6172 17 17 799 • envsales-gmbh@licor.com • envsupport-gmbh@licor.com

LI-COR Ltd., United Kingdom

Serving Denmark, Finland, Ireland, Norway, Sweden, and UK.
LI-COR Biosciences UK Ltd. • St. John's Innovation Centre • Cowley Road • Cambridge • CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422102 • Fax: +44 (0) 1223 422105 • envsales-UK@licor.com • envsupport-UK@licor.com

LI-COR Distributor Network: www.licor.com/env/distributors

© 2015 LI-COR, Inc. LI-COR is a registered trademarks of LI-COR, Inc., in the United States and other countries.
All other trademarks belong to their respective owners. For patent information, visit www.licor.com/patents