



## Soil Gas Flux Publications List

Citations related to the use of LI-COR's LI-8100/A Automated Soil Gas Flux System

*This document contains a sampling of recent publications that reference LI-COR instrumentation. 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.*

January 1, 2012 to June 30, 2015

Volume 1

- Amonette J. E., et al. (2013). **Measurement of advective soil gas flux: results of field and laboratory experiments with CO<sub>2</sub>**. *Environmental Earth Sciences* 70(4): 1717–1726.
- Anderson-Teixeira, K. J., et al. (2013). **Altered belowground carbon cycling following land-use change to perennial bioenergy crops**. *Ecosystems* 16(3): 508–520.
- Andrew, et al. (2014). **Aspen-associated mycorrhizal fungal production and respiration as a function of changing CO<sub>2</sub>, O<sub>3</sub> and climatic variables**. *Fungal Ecology* 10: 70–80.
- Anish, et al. (2014). **Organic management impacts on micro-environment in cassava**. *Journal of Root Crops* 40(1): 102–104.
- Batson, et al. (2015). **Soil greenhouse gas emissions and carbon budgeting in a short-hydroperiod floodplain wetland**. *Journal of Geophysical Research* 120(1): 77–95. January.
- Bicalho, et al. (2014). **Spatial variability structure of soil CO<sub>2</sub> emission and soil attributes in a sugarcane area**. *Agriculture, Ecosystems and Environment* 189: 206–215.
- Bortolotto, et al. (2015). **Soil carbon dioxide flux in a no-tillage winter system**. *African Journal of Agricultural Research* 10(6): 450–457.
- Bossa, A. Y., and B. Diekkruger (2014). **Spatio-temporal variability of soil respiration in a spruce-dominated headwater catchment in western Germany**. *Biogeosciences* 11: 4235–4249.
- Bowne, D. R., and E. R. Johnson (2012). **Comparison of soil carbon dioxide efflux between residential lawns and corn fields**. *Soil Science Society of American Journal* 77(3): 856–859.
- Bremenfeld, S., et al. (2013) **Effect of interrill erosion, soil crusting and soil aggregate breakdown in situ CO<sub>2</sub> effluxes**. *CATENA* 104: 14–20.
- Broeckx, L. S., et al. (2013). **The effect of a dry spring on seasonal carbon allocation and vegetation dynamics in a poplar bioenergy plantation**. *Global Change Biology: Bioenergy* 6(5): 473–487.
- Brown, K. A., et al. (2015). **Inorganic carbon system dynamics in a landfast Arctic sea ice during the early-melt period**. *Journal of Geophysical Research: Oceans* 120(5): 3542–3566.
- Bruun, E. W., et al. (2012). **Effects of slow and fast pyrolysis biochar on soil C and N turnover dynamics**. *Soil Biology and Biochemistry* 46: 73–79.
- Campoe, O. C., et al. (2012) **Stand-level patterns of carbon fluxes and partitioning in a Eucalyptus grandis plantation across a gradient of productivity, in São Paulo State, Brazil**. *Tree Physiology* 32(6): 696–706.
- Carman, C. H., et al. (2014). **Update on soil CO<sub>2</sub> Flux Monitoring at the Illinois Basin – Decatur Project, USA**. *Energy Procedia* 63: 3869–3880.
- Carvalho, J. V. D., et al. (2013). **CO<sub>2</sub> – C losses and carbon quality of selected Maritime Antarctic soils**. *Antarctic Science* 25(1): 11–18.
- Chen, S., et al. (2012). **Effects of simulated acid rain on soil CO<sub>2</sub> emissions in a secondary forest in subtropical China**. *Geoderma* 189–190: 65–71.
- Cohen, L. R., et al. (2014). **Measuring diurnal cycles of evapotranspiration in the Arctic with an automated chamber system**. *Ecohydrology* 8(4): 652–659.
- Cui, J., et al. (2012). **Changes in soil carbon sequestration and soil respiration following afforestation on paddy fields in north subtropical China**. *Journal of Plant Ecology* 6(3) 240–252.
- Daigh, A. L., et al. (2014). **Spatial and temporal dynamics of soil-surface carbon dioxide emissions in bioenergy corn rotations and reconstructed prairies**. *Soil Science Society of American Journal* 78(4): 1338–1350.

- Epron, D., et al. (2012). **Do changes in carbon allocation account for the growth response to potassium and sodium applications in tropical Eucalyptus plantations?** *Tree Physiology* 32(6): 667–679.
- Escobar, C., et al. (2015). **Biocrusts modulate warming and rainfall exclusion effects on soil respiration in a semi-arid grassland.** *Soil Biology and Biochemistry* 80: 9–17.
- Etzold, S., et al. (2013). **Long-term stem CO<sub>2</sub> concentration measurements in Norway spruce in relation to biotic and abiotic factors.** *New Phytologist* 197: 1173–1184.
- Fa, K., et al. (2015). **CO<sub>2</sub> absorption of sandy soil induced by rainfall pulses in a desert ecosystem.** *Hydrological Processes* 29(8): 2043–2051.
- Fan, J., et al. (2012). **Effects of precipitation pulses on water and carbon dioxide fluxes in two semiarid ecosystems: measurements and modeling.** *Environmental Earth Sciences* 67: 2315–2324.
- Fiener, P., et al. (2012). **Spatial variability of soil respiration in a small agricultural watershed – Are patterns of soil redistribution important?** *CATENA* 94: 3–16.
- Figueiredo, E. B., et al. (2014). **Short-term CO<sub>2</sub>-C emissions from soil prior to sugarcane (*Saccharum spp.*) replanting in southern Brazil.** *Global Change Biology: Bioenergy* 7(2): 316–327.
- Fisher, M. L., et al. (2012). **Carbon, water, and heat flux responses to experimental burning and drought in a tallgrass prairie.** *Agricultural and Forest Meteorology* 166: 169–174.
- Flores, J. P. and B. Tracy (2012). **Impacts of winter hay feeding on pasture soils and plants.** *Agriculture, Ecosystems, and Environment* 149: 30–36.
- Gabriel, C., et al. (2014). **Investigating the role of moisture as an environmental constraint in the decomposition of shallow and deep mineral soil organic matter of a temperate coniferous soil.** *Soil Biology and Biochemistry* 68: 373–384.
- Gelsomino, A., et al. (2012). **Effects of carbon sequestration methods on soil respiration and root systems in microcosm experiments and in vitro studies.** *Carbon Sequestration in Agricultural Soils*. 261–290.
- Gomez-Casanovas, N., et al. (2012). **Net ecosystems exchange modifies the relationship between the autotrophic and heterotrophic components of soil respiration with abiotic factors in prairie grasslands.** *Global Change Biology* 18(8): 2532–2545.
- Gomez-Casanovas, N., et al. (2013). **Gap filling strategies and error in estimating annual soil respiration.** *Global Change Biology* 19(6): 1941–1952.
- Gong, Y. M., et al. (2014). **Response of carbon dioxide emissions to sheep grazing and N application in an alpine grassland – Part 1: Effect of sheep grazing.** *Biogeosciences* 11: 1743–1750.
- Gonzalez, S., et al. (2014). **Climate factors mediate soil respiration dynamics in Mediterranean agricultural environments: an empirical approach.** *Soil Research* 52(6): 543–553.
- Graf, A., et al. (2012). **Analyzing spatiotemporal variability of heterotrophic soil respiration at the field scale using orthogonal functions.** *Geoderma* 181: 91–101.
- Groh, T. A., et al. (2014). **Nitrogen removal and greenhouse gas emissions from constructed wetlands receiving tile drainage water.** *Journal of Environmental Quality* 44(3): 1001–1010.
- Grossiord, C., et al. (2012). **Transpiration alters the contribution of autotrophic and heterotrophic components of soil CO<sub>2</sub> efflux.** *New Phytologist* 194(3): 647–653.
- Hagedorn, F., et al. (2012). **Nitrogen addition alters mineralization dynamics of <sup>13</sup>C-depleted leaf and twig litter and reduces leaching of older DOC from mineral soil.** *Global Change Biology* 18(4): 1412–1427.
- Hagedorn, F. and O. Joos (2014). **Experimental summer drought reduces soil CO<sub>2</sub> effluxes and DOC leaching in Swiss grassland soils along an elevational gradient.** *Biogeochemistry* 117: 395–412.
- Hamerlynck, E. P., et al. (2013). **Consequences of cool-season drought-induced plant mortality to Chihuahuan desert grassland ecosystems and soil respiration dynamics.** *Ecosystems* 16: 1178–1191.
- Han, Y., et al. (2012). **Effects of mowing and nitrogen addition on soil respiration in three patches in an oldfield grassland in Inner Mongolia.** *Journal of Plant Ecology* 5(2): 219–228.

- Han, W., et al. (2014). **Accounting for time-variable soil porosity improves the accuracy of the gradient method for estimating soil carbon dioxide production.** *Soil Science Society of America Journal* 78(4): 1426–1433.
- Heinemeyer, A., et al. (2012) **Partitioning of soil CO<sub>2</sub> flux components in a temperate grassland ecosystem.** *European Journal of Soil Science* 63(2): 249–260.
- Hu, Z., et al. (2013). **Soil respiration and N<sub>2</sub>O flux response to UV-B radiation and straw incorporated in a soybean-winter wheat rotation system.** *Water Air Soil Pollution* 224: 1394.
- Hu, Y., et al. (2014). **Variability in soil microbial community and activity between coastal and riparian wetlands in the Yangtze River estuary – potential impacts on carbon sequestration.** *Soil Biology and Biochemistry* 70: 221–228.
- Hu, Z., et al. (2014). **Effects of harvest residue management on soil carbon and nitrogen processes in a Chinese fir plantation.** *Forest Ecology and Management* 326: 163–170.
- Ibrahim, M., et al. (2015). **Changes of CO<sub>2</sub> emissions and labile organic carbon as influenced by rice straw and different water regimes.** *International Journal of Environmental Science and Technology* 12: 263–274.
- Jensen, A. E., et al. (2014). **Variations in soil carbon dioxide efflux across a thaw slump chronosequence in northwestern Alaska.** *Environmental Research Letters* 9: 025001.
- Jiang, J., et al. (2015). **Changes in temperature sensitivity of soil respiration in the phases of a three-year crop rotation system.** *Soil and Tillage Research* 150: 139–146.
- Jiang, M., et al. (2015). **Improved understanding of climate change impact to Pennsylvania dairy pastures.** *Crop Science* 55(2): 934–949.
- Jones, D. G., et al. (2014). **Monitoring of near surface gas seepage from a shallow injection experiment at the CO<sub>2</sub> field lab, Norway.** *International Journal of Greenhouse Gas Control* 28: 300–317.
- Kaplan, H., et al. (2014). **Impact of trace metal contamination and in situ remediation in microbial diversity and respiratory activity of heavily polluted kastanozems.** *Biology and Fertility of Soils* 50(5): 735–744.
- Keidel, L., et al. (2014). **Positive feedback of elevated CO<sub>2</sub> on soil respiration in late autumn and winter.** *Biogeosciences Discussions* 11: 8749–8787.
- Kurth, V. J., et al. (2014). **Initial soil respiration response to biomass harvesting and green-tree retention in aspen-dominated forests of the Great Lakes region.** *Forest Ecology and Management*. 15: 342–352.
- Lay, C., et al. (2012). **Microbial diversity and activity in hypersaline high Arctic spring channels.** *Extremophiles* 16: 177–191.
- Lewicki, J. L., et al. (2014). **Multi-scale observations of the variability of magnetic CO<sub>2</sub> emissions, Mammoth Mountain, CA, USA.** *Journal of Volcanology and Geothermal Research* 284: 1–15.
- Li, D., et al. (2013). **Contrasting response of heterotrophic and autotrophic respiration to experimental warming in a winter annual-dominated prairie.** *Global Change Biology* 19: 3553–3564.
- Liu, D., et al. (2013). **Effect of rain enrichment on soil respiration of *Nitraria sphaerocarpa* community in a hyperarid area.** *Sciences in Cold and Arid Regions* 5(6): 722–732.
- Liu, Y., et al. (2014). **Variation in soil respiration under the tree canopy in a temperate mixed forest, central China, under different soil water conditions.** *Ecological Research* 29: 133–142.
- Loisy, C., et al. (2013). **The CO<sub>2</sub>-Vadose Project: Dynamics of the natural CO<sub>2</sub> in a carbonate vadose zone.** *International Journal of Greenhouse Gas Control*. 14: 97–112.
- Ma, J., et al. (2012). **The response of CO<sub>2</sub> flux to rain pulses at a saline desert.** *Hydrological Processes* 26(26): 4029–4037.
- Ma, Y., et al. (2013). **Effect of clear-cutting silviculture on soil respiration in a subtropical forest of China.** *Journal of Plant Ecology* 6(5): 335–348.
- Ma, Y., et al. (2014). **Stand ages regulate the response of soil respiration to temperature in a *Larix principis-rupprechtii* plantation.** *Agriculture and Forest Meteorology*. 184: 179–187.
- Maestre, F. T., et al. (2013). **Changes in biocrust cover drive carbon cycle responses to climate change in drylands.** *Global Change Biology* 19(12): 3835–3847.

- Marasco, D. E., et al. (2014). **Quantifying evapotranspiration from urban green roofs: A comparison of chamber measurements with commonly used predictive methods.** *Environmental Science and Technology*. 48(17): 10273–10281.
- Martin, J. G., et al. (2012). **High-frequency analysis of the complex linkage between soil CO<sub>2</sub> fluxes, photosynthesis and environmental variables.** *Tree Physiology* 32(1): 49.
- Mathiba, M. and Awuah-Offei, K. (2015). **Spatial autocorrelation of soil CO<sub>2</sub> fluxes on reclaimed mine lands.** *Environmental Earth Sciences* 73(12): 8287–8297.
- Mathiba, M., et al. (2015). **Influence of elevation, soil temperature and soil moisture content on reclaimed mine land soil CO<sub>2</sub> fluxes.** *Environmental Earth Sciences*. 73(10): 6131–6143.
- Matías, L., et al. (2012). **Effect of simulated climate change on soil respiration in a Mediterranean-type ecosystem: rainfall and habitat type are more important than temperature or the soil carbon pool.** *Ecosystems* 15: 299–310.
- Mazot, A., et al. (2015). **CO<sub>2</sub> Degassing from Volcanic Lakes.** *Volcanic Lakes (Advances in Volcanology)*. 341–354.
- Miao, G., et al. (2013). **The effect of water table fluctuation on soil respiration in a lower coastal plain forested wetland in the southeastern U.S.** *Journal of Geophysical Research: Biogeosciences* 118 (4). 1748–1762.
- Mukhopadhyay, S., and S. K. Maiti (2012). **Soil CO<sub>2</sub> Flux in grassland, afforested land and reclaimed coalmine overburden dumps: A case study.** *Land Degradation and Development*. 25(3): 216–227.
- Nagano, T., et al. (2013). **Subsidence and soil CO<sub>2</sub> efflux in tropical peatland in southern Thailand under Various water table and management conditions.** *Mires and Peat* 11: 1–20.
- Noormets, A., et al. (2012). **The role of harvest residue in rotation cycle carbon in loblolly pine plantations. Respiration partitioning approach.** *Global Change Biology* 18(10): 3186–3201.
- Nottingham, A. T., et al. (2012). **Soil properties in tropical montane cloud forest influence estimates of soil CO<sub>2</sub> efflux.** *Agricultural and Forest Meteorology* 166: 215–220.
- Pacaldo, R. S., et al. (2014). **Soil CO<sub>2</sub> effluxes, temporal and spatial variations, and root respiration in shrub willow biomass crops fields along a 19-year chronosequence as affected by regrowth and removal treatments.** *Global Change Biology: Bioenergy* 6(5): 488–498.
- Pacaldo, R. S., et al. (2014). **Carbon sequestration in fine roots and foliage offsets soil CO<sub>2</sub> effluxes along a 19-year chronosequence of shrub willow biomass crops.** *Bioenergy Resources* 7: 769–776.
- Pang, Z., et al. (2012). **Environmental relationships with the interannual and seasonal variation of soil respiration in a cedar (*Platycladus orientalis*) plantation in northern China.** *Canadian Journal of Forest Research* 42(5): 934–942.
- Parks, M. M., et al. (2013). **Distinguishing contributions to diffuse CO<sub>2</sub> emissions in volcanic areas from magmatic degassing and thermal decarbonation using soil gas 222Rn- $\delta^{13}C$  systematics: Application to Santorini volcano, Greece.** *Earth and Planetary Science Letters* 377: 180–190.
- Perez-Quezada, J. F., et al. (2012). **Effects of afforestation on soil respiration in an arid shrubland in Chile.** *Journal of Arid Environments* 83: 45–53.
- Potter, C. (2012). **Net primary production and carbon cycling in coast redwood forests of central California.** *Open Journal of Ecology* 2(3): 147–153.
- Pronost, J., et al. (2015). **CO<sub>2</sub>-depleted warm air venting from chrysotile milling waste (Thetford Mines, Canada): Evidence for in-situ carbon capture from the atmosphere.** *Geology* 40(3): 275–278.
- Quin, S., et al. (2014). **Restoration of upland heath from a graminoid- to a *Calluna vulgaris*-dominated community provides a carbon benefit.** *Agriculture, Ecosystems and Environment* 185: 133–143.
- Rezanezhad, F., et al. (2014). **Water table fluctuations and soil biogeochemistry: An experimental approach using an automated soil column system.** *Journal of Hydrology* 509: 245–256.
- Richards, A. E., et al. (2012). **Fire interacts with season to influence soil respiration in tropical savannas.** *Soil Biology and Biochemistry* 53: 90–98.

- Riederer, M., et al. (2014). **Net ecosystem CO<sub>2</sub> exchange measurements by the closed chamber method and the eddy covariance technique and their dependence on atmospheric conditions.** *Atmospheric Measurement Techniques* 7: 1057–1064.
- Roland, M., et al. (2015). **Importance of nondiffusive transport for soil CO<sub>2</sub> efflux in a temperate mountain grassland.** *Journal of Geophysical Research: Biogeosciences* 120(3): 502–512.
- Ruehr, N. K., et al. (2012). **Effects of water availability on carbon and water exchange in a young ponderosa pine forest: Above- and belowground responses.** *Agricultural and Forest Meteorology* 164: 136–148.
- Ruehr, N. K., et al. (2014). **Effects of heat and drought on carbon and water dynamics in a regenerating semi-arid pine forest: a combined experimental and modeling approach.** *Biogeosciences* 11: 4139–4156.
- Russell, A. E. (2014). **Unexpected effects of chitin, cellulose, and lignin addition on soil dynamics in a wet tropical forest.** *Ecosystems* 17: 918–930.
- Rutledge, S., et al. (2014). **CO<sub>2</sub> emissions following cultivation of a temperate permanent pasture.** *Agriculture, Ecosystems, and Environment* 184: 21–33.
- Ryals, R. and Silver, W. L. (2013). **Effects of organic matter amendments on net primary productivity and greenhouse gas emissions in annual grasslands.** *Ecological Applications* 23: 46–59.
- Santos, E., et al. (2014). **Temporal dynamics of oxygen isotope compositions of soil and canopy CO<sub>2</sub> fluxes in a temperate deciduous forest.** *Journal of Geophysical Research: Biogeosciences* 119(5): 996–1013.
- Satakhum, D., et al. (2013). **Soil CO<sub>2</sub> efflux and soil carbon balance of a tropical rubber plantation.** *Ecological Resources* 28: 969–979.
- Schatz, J. D., et al. (2012). **Spatial patterns of soil surface C flux in experimental canopy gaps.** *Ecosystems* 15: 616–623.
- Shao, R., et al. (2014). **Nitrogen fertilization increase soil carbon dioxide efflux of winter wheat field: A case study in Northwestern China.** *Soil and Tillage Research* 143: 164–171.
- Silva-Olaya, A. M., et al. (2013). **Carbon dioxide emissions under different soil tillage systems in mechanically harvested sugarcane.** *Environmental Research Letters* 8(1).
- Singh, K., et al. (2012). **Ecological restoration of degraded sodic lands through afforestation and cropping.** *Ecological Engineering* 43: 70–80.
- Singh, K., et al. (2012). **Changes in physio-chemical, microbial and enzymatic activities during restoration of degraded sodic land: Ecological suitability of mixed forest over monoculture plantation.** *CATENA* 96: 57–67.
- Smith, S. W., et al. (2014). **Root traits predict decomposition across a landscape-scale grazing experiment.** *New Phytologist* 203: 851–862.
- Song, W., et al. (2012). **Vegetation cover and rain timing co-regulate the response of soil CO<sub>2</sub> efflux to rain increase in an arid desert ecosystem.** *Soil Biology and Biochemistry* 49: 114–123.
- Song, W., et al. (2015). **Simulated rain addition modifies diurnal patterns and temperature sensitivities of autotrophic and heterotrophic soil respiration in an arid desert ecosystem.** *Soil Biology and Biochemistry* 82: 143–152.
- Soong, J. L. and M. F. Cotrufo (2015). **Annual burning of a tallgrass prairie inhibits C and N cycling in soil, increasing recalcitrant pyrogenic organic matter storage while reducing N availability.** *Global Change Biology*. 21: 2321–2333.
- Sun, Z., et al. (2014). **The effect of nitrogen addition on soil respiration from a nitrogen-limited forest soil.** *Agricultural and Forest Meteorology* 197: 103–110.
- Thomazini, A., et al. (2014). **Spatial Variability of CO<sub>2</sub> Emissions from Newly Exposed Paraglacial Soils at a Glacier Retreat Zone on King George Island, Maritime Antarctica.** *Permafrost and Periglacial Processes* 25(4): 233–242.
- Troxler, T. J., et al. (2015). **Component-specific dynamics of riverine mangrove CO<sub>2</sub> efflux in the Florida coastal Everglades.** *Agricultural and Forest Meteorology*. Available online: February, 2015.
- Versini, A., et al. (2013). **The manipulation of organic residues affects tree growth and heterotrophic CO<sub>2</sub> efflux in a tropical *Eucalyptus* plantation.** *Forest Ecology and Management* 301: 79–88.

- Wall, G. W., et al. (2013). **Infrared warming affects intrarow soil carbon dioxide efflux during vegetative growth of spring wheat.** *Agronomy Journal* 105(3): 607–618.
- Wang, J. (2012). **Annual soil CO<sub>2</sub> efflux in a wet meadow during active layer freeze-thaw changes on the Qinghai-Tibet Plateau.** *Environmental Earth Science* 69(3) 855–862.
- Wang, C. (2013). **Seasonality of soil CO<sub>2</sub> efflux in a temperate forest: Biophysical effects of snowpack and spring freeze-thaw cycles.** *Agriculture and Forest Meteorology* 177: 83–92.
- Wang, W., et al. (2014). **Modeling the contribution of abiotic exchange to CO<sub>2</sub> flux in alkaline soils of arid areas.** *Journal of Arid Land* 6(1): 27–36.
- Wang, Y., et al. (2014). **Non-additive effects of litter-mixing on soil carbon dioxide efflux from poplar-based agroforestry systems in the warm temperate region of China.** *Agroforestry Systems* 88(2): 193–203.
- Wang, Y., et al. (2014). **Non-growing-season soil respiration is controlled by freezing and thawing processes in the summer monsoon-dominated Tibetan alpine grassland.** *Global Biogeochemical Cycles* 28(10): 1081–1095.
- Wang, J., et al. (2014). **Variations in leaf litter decomposition across contrasting forest stands and controlling factors at local scale.** *Journal of Plant Ecology*. Advance access: October, 2014.
- Warren, C. R., (2014). **Response of osmolytes in soil to drying and rewetting.** *Soil Biology and Biochemistry* 70: 22–32.
- Wei, S., et al. (2014). **Effect of soil temperature and soil moisture on CO<sub>2</sub> flux from eroded landscape positions on black soil in Northeastern China.** *Soil and Tillage Research* 144: 119–125.
- Xia, J., et al. (2012). **The effects of warming-shifted plant phenology on ecosystem carbon exchange are regulated by precipitation in a semi-arid grassland.** *PLoS one* 7(2): e32088.
- Xiao, C., et al. (2007). **Irrigation and enhanced soil carbon input effects on below-ground carbon cycling in semiarid temperate grasslands.** *New Phytologist* 174: 835–846.
- Xie, t., P. Su (2012). **Canopy and leaf photosynthetic characteristics and water use efficiency of sweet sorghum under drought stress.** *Russian Journal of Plant Physiology*. 59(2): 224–234.
- Xu, X., et al. (2015). **Plant Community structure regulates responses of prairie soil respiration to decadal experimental warming.** *Global Change Biology*. First published online: June, 2015.
- Zanchi, F. B., et al. (2014). **Soil CO<sub>2</sub> exchange in seven pristine Amazonian rain forest sites in relation to soil temperature.** *Agricultural and Forest Meteorology* 192: 96–107.
- Zanotelli, D., et al. (2013). **Net Primary productivity, allocation pattern and carbon use efficiency in an apple orchard assessed by integrating eddy covariance, biometric and continuous soil chamber measurements.** *Biogeosciences* 10: 3089–3108.
- Zhang, Y., et al. (2012). **Chemistry of decomposing mulching materials and the effect on soil carbon dynamics under *Phyllostachys praecox* bamboo stand.** *Journal of Soils and Sediments* 13: 24–33.
- Zhang, L., et al. (2012). **Soil carbon dioxide flux from shelterbelts in farmland in temperate arid region, northwest China.** *European Journal of Soil Biology* 48: 24–31.
- Zhang, X., et al. (2012). **Effects of raised-bed planting for enhanced summer maize yield on rhizosphere soil microbial functional groups and enzyme activity in Henan Province, China.** *Field Crops Research* 130: 28–37.
- Zhang, T., et al. (2013). **Responses of seasonal and diurnal soil CO<sub>2</sub> effluxes to land-use change from paddy fields to Lei bamboo (*Phyllostachys praecox*) stands.** *Atmospheric Environment*. 77: 856–864.
- Zheng, X., et al. (2014). **Soil CO<sub>2</sub> efflux along an elevation gradient in Qinghai spruce forests in the upper reaches of the Heihe River, northwest China.** *Environmental Earth Sciences* 71(5): 2065–2076.
- Zhou, Z., et al. (2013). **Predicting soil respiration using carbon stock in roots, litter and soil organic matter in forests of loess Plateau in China.** *Soil Biology and Biochemistry* 57: 135–143.
- Zimmermann, M. and M.I. Bird (2012). **Temperature sensitivity of tropical forest soil respiration increase along an altitudinal gradient with ongoing decomposition.** *Geoderma* 187: 8–15.
- Zimmermann, M., et al. (2012). **Rapid degradation of pyrogenic carbon.** *Global Change Biology* 18: 3306–3316.

Please contact us with any corrections or potential additions to this list.

Tell us about your research by visiting [www.licor.com/case-study](http://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](mailto:envsales@licor.com) • [envsupport@licor.com](mailto:envsupport@licor.com) • [www.licor.com/env](http://www.licor.com/env)

### ***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](mailto:envsales-gmbh@licor.com) • [envsupport-gmbh@licor.com](mailto: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](mailto:envsales-UK@licor.com) • [envsupport-UK@licor.com](mailto:envsupport-UK@licor.com)

**LI-COR Distributor Network:** [www.licor.com/env/distributors](http://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](http://www.licor.com/patents)

The LI-COR logo is displayed in a bold, italicized, black sans-serif font. The letters 'LI' are connected to 'COR', and there is a registered trademark symbol (®) at the end.