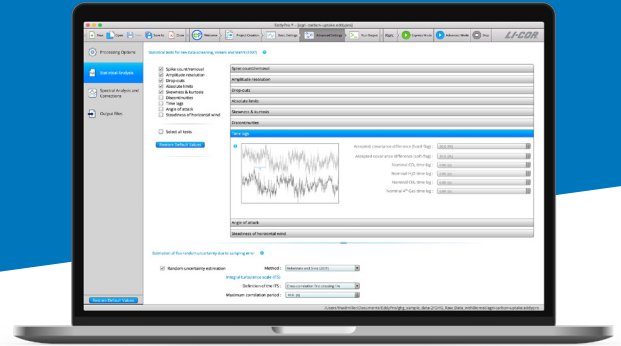




EddyPro[®] Software

The Community Standard



EddyPro Software is the only flux processing software developed in collaboration with FLUXNET, AmeriFlux, ICOS, and the eddy covariance research community. EddyPro uses the most accepted and cited algorithms to compute fully-processed fluxes of water vapor, carbon dioxide, methane, other trace gases, and energy, using the eddy covariance method.

With over 7,000 downloads in 176 countries, EddyPro has been widely adopted by eddy covariance researchers and networks as their standard processing software. Version 7 includes new variables, algorithms, and outputs not provided by any other software. Achieve fully processed fluxes on site and in real time with EddyPro running on the SmartFlux[®] System.

To learn more about EddyPro and its features visit licor.com/eddypro.

Key Features

- Outputs conform with world standards in variable naming, formatting, and unit structure for easy comparison and network submission
- Fully processed fluxes on site and in real time with EddyPro running on the SmartFlux System
- EddyPro version 7 includes variables, algorithms, and outputs only offered by LI-COR
- EddyPro is freely available and open source
- User-friendly interface, updated with the latest data processing methods, supported and documented by LI-COR

EddyPro® Software

System Requirements

EddyPro is compatible with computers running the Windows® and MacOS® operating system.

Foundational References

- Burba et al. (2008) – Global Change Biology, 14:1854–1876.
- Burba et al. (2012) – Global Change Biology, 18:385-399.
- Finkelstein and Sims (2001) – Journal of Geophysical Research, 106:3,503 – 3,509.
- Foken (2003) – Angewandte Meteorologie, Mikrometeorologische Methoden, 289 pp.
- Foken et al. (2004) – Handbook of micrometeorology: A guide for surface flux measurements, 81-108.
- Fratini et al. (2012) – Agricultural and Forest Meteorology, 165: 53-63
- Göckede et al. (2004) – Agricultural and Forest Meteorology, 127: 175-188.
- Hsieh et al. (2000) – Advances in Water Resources, 23: 765-772.
- Horst (1997) – Boundary Layer Meteorology, 82: 219-233.
- Horst and Lenschow (2009) – Boundary Layer Meteorology, 130: 275-300.
- Ibrom et al. (2007a) – Tellus Series B-Chemical and Physical Meteorology, 59:937-946.
- Ibrom et al. (2007b) – Agricultural and Forest Meteorology, 147: 140-156.
- Kljun et al. (2004) – Boundary Layer Meteorology, 112: 503-523.
- Kormann and Meixner (2001) – Boundary Layer Meteorology, 99:207–224.
- Mann and Lenschow (1994) – Journal of Geophysical Research, 99:14,519–14,526.
- Mauder and Foken (2004) - Documentation and Instruction Manual of the Eddy Covariance Software Package TK2, Arbeitsergebnisse Nr. 26.
- McDermitt et al. (2011) – Applied Physics B: Laser and Optics, 102: 391-405.
- Moncrieff et al. (1997) – Journal of Hydrology, 188-189: 589-611.
- Moncrieff et al. (2004) – Handbook of micrometeorology, 7-31.
- Nakai et al. (2006) – Agricultural and Forest Meteorology, 136: 19-30.
- Nakai and Shimoyama (2012) – Agricultural and Forest Meteorology, 162-163: 14-26
- Van Dijk et al. (2004) – Meteorology and Air Quality Group, Wageningen, the Netherlands, 99 pp.
- Vickers and Mahrt (1997) – Journal of Atmospheric and Oceanic Technology, 14: 512-526.
- Webb et al. (1980) – Quarterly Journal of the Royal Meteorological Society, 106: 85-100.
- Wilczak et al. (2001) - Boundary Layer Meteorology, 99: 127-150.