

Systems Integration

Bringing It All Together

Transform your eddy covariance site into an entire ecosystem research site. LI-COR eddy covariance systems complement or integrate additional biological flux data from other instruments and components to provide a more complete picture of the activity in your ecosystem.



Soil Flux Data

The LI-8100A Automated Soil CO₂ Flux system with LI-8150 multiplexer and long-term chambers is ideal for assessing the temporal and spatial variability of soil CO₂ fluxes at your eddy covariance site. Including soil flux data with your eddy covariance data can provide a more comprehensive understanding of ecosystem behavior. Soil data from the LI-COR Soil Flux System can be integrated and time synchronized using EddyPro[®] Software running on the SmartFlux[®] System for real-time processing with your eddy covariance flux data.

Learn more about the LI-8100A at licor.com/soil.



Photosynthesis Data

Including photosynthetic data can help you build your carbon budget with consideration of carbon sequestration and gas fluxes from plants. The LI-6800 Portable Photosynthesis System quickly delivers superior photosynthetic gas exchange and chlorophyll fluorescence measurements.

See the latest advancements in photosynthesis research at licor.com/6800.



Leaf Area Index

Add details of your ecosystem's dynamic canopy with leaf area index measurements. Leaf Area Index (LAI) describes an ecosystem's dynamic canopy and is often required when submitting eddy covariance data to worldwide networks. It plays an important role in carbon, water, and energy fluxes. The LAI-2200C is a standalone instrument for manual, non-destructive LAI measurements.

For more information about the LAI-2200C visit licor.com/2200C.



Data Site Images

For daily images of your site, the PhenoCam is a robust outdoor camera that integrates with LI-COR eddy covariance systems. You can view PhenoCam images on FluxSuite® Software or download them from your site.

For details on adding a PhenoCam to your eddy covariance site visit licor.com/biomet.