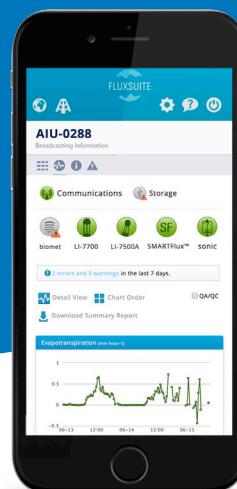


# FluxSuite® Software

## Always Online Research Site Monitoring



You cannot be everywhere at once, but your eddy covariance site can. FluxSuite® Software lets you put your eddy covariance sites online so you can monitor fluxes, biomet data, and status information from anywhere, at any time, on any device, whether that's a smartphone, tablet or computer.

FluxSuite is a secure web service that provides status information and on-site, real-time results, processed by EddyPro® running on the SmartFlux® System, from your eddy covariance site. Receive email notifications any time an instrument triggers an alert, so you can go to the site prepared to solve the problem.

Add as many stations as you want. View data from multiple sensors at any station and assign collaborators to simplify site management.

To learn more about FluxSuite and its features visit [fluxsuite.com](http://fluxsuite.com).

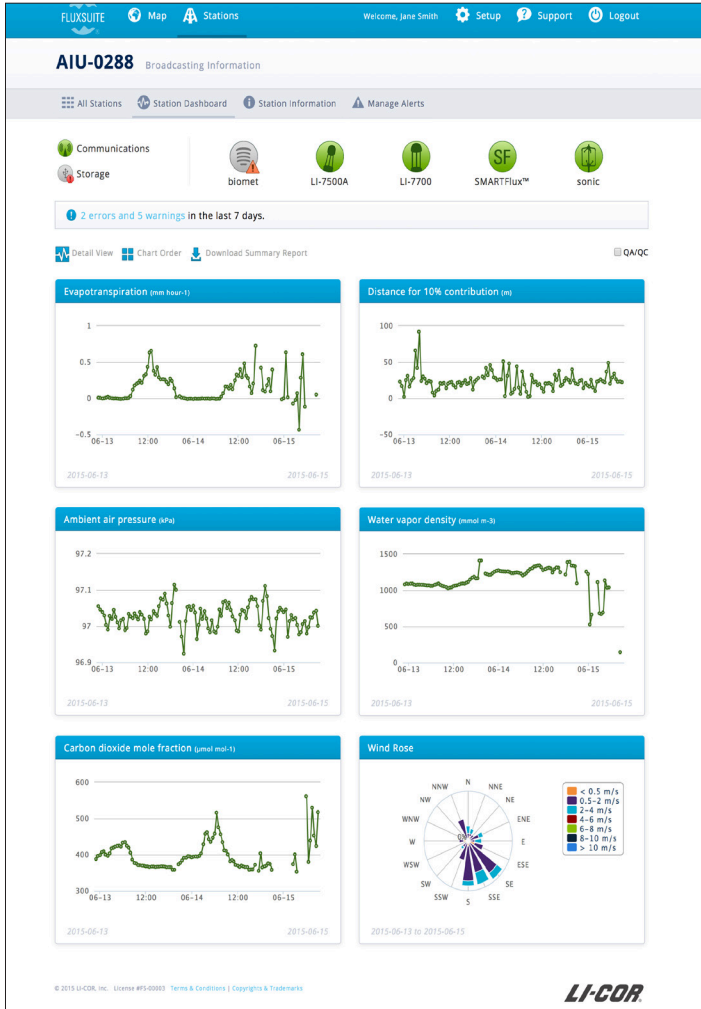
### Key Features

- Monitor sensor performance and condition to maximize data collection
- Share real-time results with team members or collaborators
- Real-time flux quality indicators
- Visit only as needed and avoid unnecessary site visits
- Reduce data loss with immediate email alerts when something needs attention
- Download and view real-time final fluxes on your smartphone, tablet, or computer

# FluxSuite® Software

## Add Convenience and Comfort to Your Research

FluxSuite Software lets you access your eddy covariance stations from anywhere in the world. Check the status of instruments, transfer raw data from your site, and configure instruments, all from your office or home. Limit the time you spend on field visits while ensuring your site continuously records data.



Station Name	Status	Abbreviation	Location	Ecosystem	Action
AIU-447	●	LERS1	LI-COR, Lincoln NE	Grassland	ⓘ
Pyrenees Conifer	●	PC	Andorra	Forest	ⓘ
Filler Station	●	FS	Titusville, FL	Scrub Oak	ⓘ
Goose Lake	●	Goose	Marshall, WI	Forest	ⓘ
AIU-0288	●	Open Path	LI-COR, Lincoln NE	Grassland	ⓘ
Serengeti	●	SER	Tanzania	Savanna	ⓘ
Zhuzhou Paddy	●	ZP	Hunan Province, China	Rice	ⓘ
Lewis and Clark	●	LC	Yankton, SD	River	ⓘ
Mobile EC	●	MEC	LERS West	Grassland	ⓘ
Sonoran	●	SD	Arizona	Desert	ⓘ

