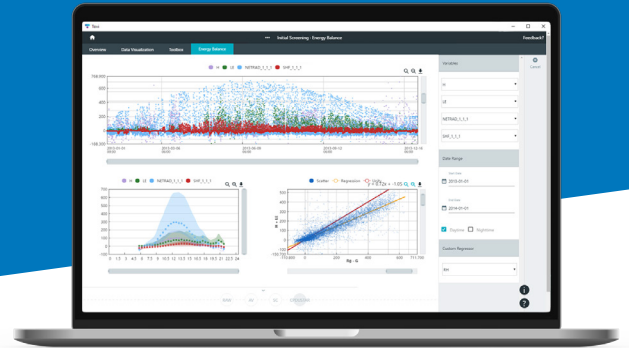




# Tovi™ Software

## Inspired Tools for Ecosystem Data Analysis



Eddy covariance data analysis does not end with fully-processed fluxes from EddyPro® Software. Tovi Software provides all the tools you need for a fully documented, traceable workflow that improves reproducibility of your research. Detailed graphs are automatically generated, and energy balance and carbon cycle dashboards help you visualize your energy budget and carbon accumulation in the ecosystem.

Tovi performs complete, scientifically-sound analysis of your EddyPro outputs and provides documentation of your processing steps for publication and collaboration in a fraction of the time it takes with other applications. Tovi integrates and processes all data to deliver reliable ecosystem insights and experimental evaluations quickly.

With Tovi, all scientific codes and algorithms used by major eddy covariance networks are combined in a single application. Full credit is given to the authoring scientists in the citation builder, which compiles a complete list of formatted citations. Tovi is continuously updated, with more tools added regularly.

To learn more about Tovi and its features visit [licor.com/tovi](https://licor.com/tovi).

### Key Features

- Quality control screening, flagging, and filtering in a single software package
- Automated weather station data retrieval
- See where your fluxes are coming from and view fluxes from different areas around your tower
- Improve data continuity by gap filling missing weather and flux data
- Flag and remove periods with unreliable flux due to low turbulence
- View only the fluxes related to your research by separating flux data into key components

# Tovi™ Software

## System Requirements

Tovi will run on most Windows® and Mac® computers. For the best performance, follow the guidelines below.

### Operating System:

64-bit Windows 7 or 64-bit Windows 10  
OSX El Capitan or newer (MacOS)

### Minimum Hardware:

Intel Core 2 - 2.0 GHz  
8 GB RAM  
1680 × 1050 pixels

### Recommended Hardware:

Intel i3 - 2.6 GHz or better  
16 GB RAM  
1920 × 1080 pixels

## Foundational References

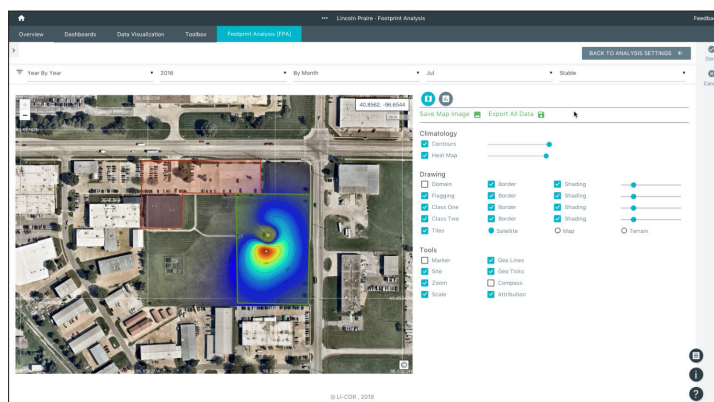
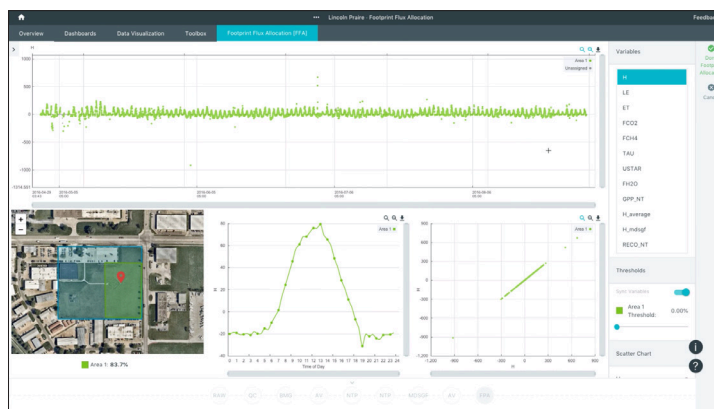
Barr, A., Richardson, A., Hollinger, D., Papale, D., Arain, M., Black, T., Bohrer, G., Dragoni, D., Fischer, M., Gu, L., Law, B., Margolis, H., McCaughey, H., Munger, W., Oechel, W. and Schaeffer, K.: Use of change-point detection for friction–velocity threshold evaluation in eddy-covariance studies, *Agricultural and Forest Meteorology*, 171, 31–45, 2013.

Charuchittipan, D., Babel, W., Mauder, M., Leps, J.-P. and Foken, T.: Extension of the averaging time in eddy-covariance measurements and its effect on the energy balance closure, *Boundary-layer meteorology*, 152(3), 303–327, doi: 10.1007/s10546-014-9922-6, 2014.

De Roo, F., Zhang, S., Huq, S. and Mauder, M.: A semi-empirical model of the energy balance closure in the surface layer, *PLoS one*, 13(12), e0209022, 2018.

Isaac, P., Cleverly, J., McHugh, I., van Gorsel, E., Ewenz, C. and Beringer, J.: OzFlux data: Network integration from collection to curation, *Biogeosciences*, 14(12), 2903–2928, doi: 10.5194/bg-14-2903-2017, 2017.

Kljun, N., Calanca, P., Rotach, M. and Schmid, H.: A simple two-dimensional parameterisation for Flux Footprint Prediction (FFP), *Geoscientific Model Development*, 8(11), 3695–3713, doi:10.5194/gmd-8-3695-2015, 2015.



Mauder, M., Cuntz, M., Drüe, C., Graf, A., Rebmann, C., Schmid, H. P., Schmidt, M. and Steinbrecher, R.: A strategy for quality and uncertainty assessment of long-term eddy-covariance measurements, *Agricultural and Forest Meteorology*, 169, 122–135, doi: <https://doi.org/10.1016/j.agrformet.2012.09.006>, 2013.

Reichstein, M., Falge, E., Baldocchi, D., Papale, D., Aubinet, M., Berbigier, P., Bernhofer, C., Buchmann, N., Gilmanov, T., Granier, A., Grünwald, T., Havránková, K., Ilvesniemi, H., Janous, D., Knohl, A., Laurila, T., Lohila, A., Loustau, D., Matteucci, G., Meyers, T., Miglietta, F., Ourcival, J. M., Pump-anen, J., Rambal, S., Rotenberg, E., Sanz, M., Tenhunen, J., Seufert, G., Vaccari, F., Vesala, T., Yakir, D. and Valentini, R.: On the separation of net ecosystem exchange into assimilation and ecosystem respiration: review and improved algorithm, *Global Change Biology*, 11(9), 1424–1439, doi:10.1111/j.1365-2486.2005.001002.x, 2005.

Seo, T.: A Microclimatological Study of Thermal Exchange at the Earth's Surface, 1958.