

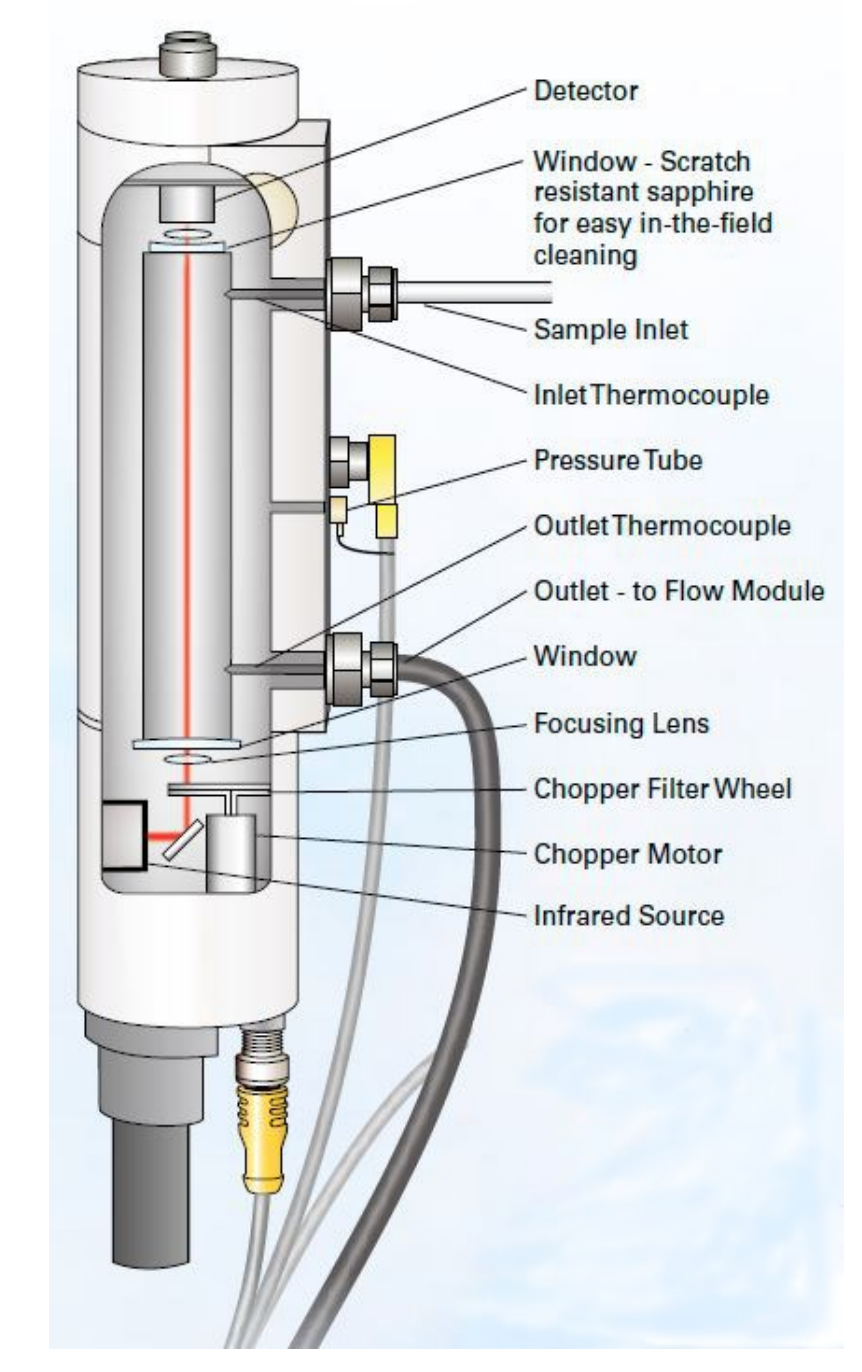
INVESTIGATION ON THE IMPORTANCE OF FAST AIR TEMPERATURE MEASUREMENTS IN THE SAMPLING CELL OF SHORT-TUBE CLOSED-PATH GAS ANALYZER FOR EDDY-COVARIANCE FLUXES

James Kathilankal, Gerardo Fratini and George Burba
LI-COR Biosciences, Lincoln, NE, USA

INTRODUCTION

- Enclosed gas analyzer design allows eddy covariance flux measurements using short intake tubes, $\leq 1\text{m}$ [1]
- Tubes with length-to-diameter ratio $< 1000:1$ do not eliminate 100% of fast temperature fluctuations [2]
- Fast temperature fluctuations affect gas density measurements and conversion to dry mole fraction [3], thus affecting flux calculations
- Computing fluxes using short-tube instruments requires high-speed temperature measurements of the air stream inside the sampling cell
- Here we examine importance of such measurements, and show flux errors resulted from not using fast temperature of air sample measured inside the cell

METHOD



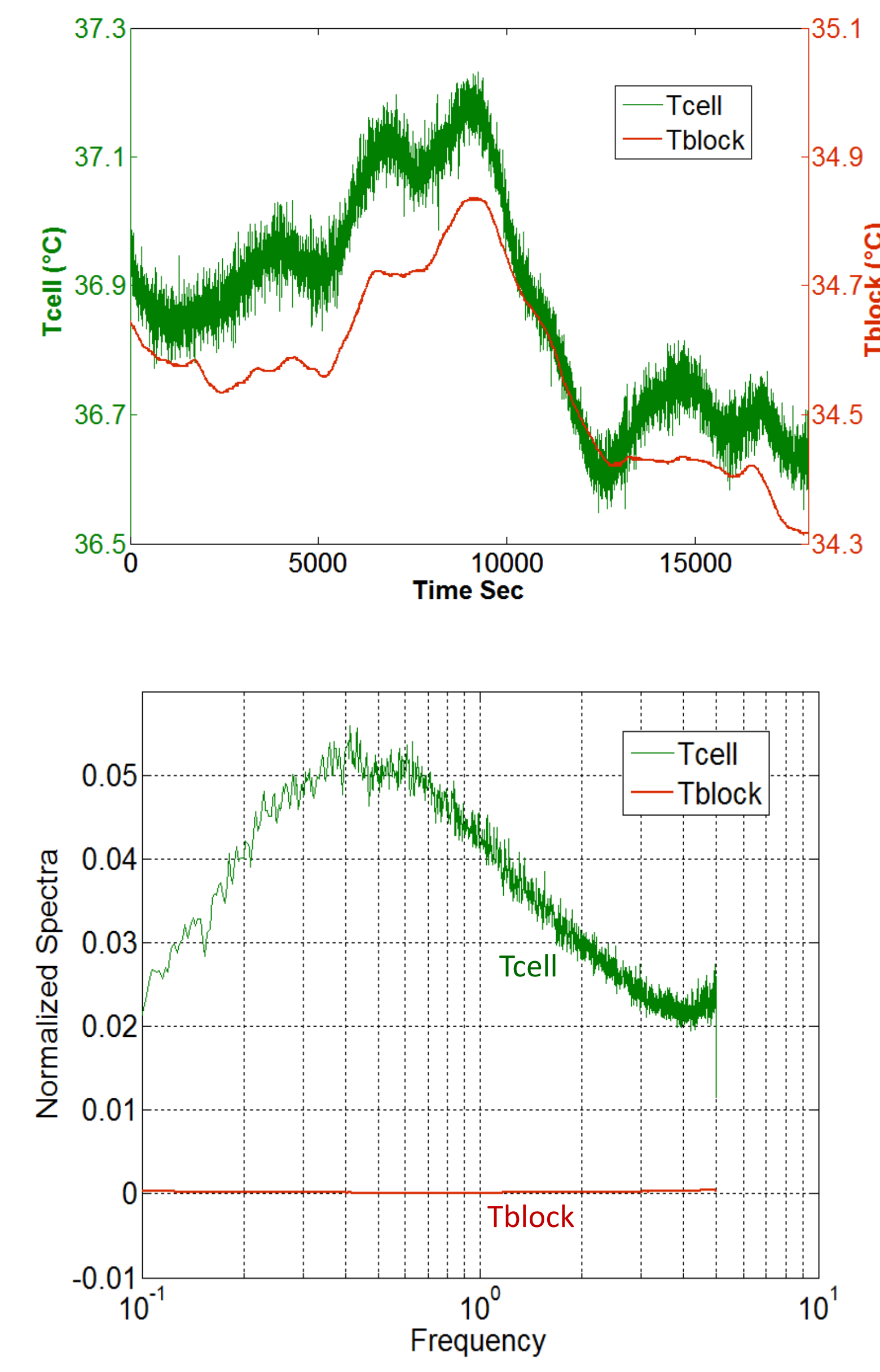
- Three LI-7200 enclosed gas analyzers with 1 m long 0.005 m diameter intake tubes were used at three experimental sites
- Each analyzer is equipped with two high-speed thermocouples weighed to output correct cell air temperature [1]
- Each analyzer also outputs slow block temperature measured in the metal body next to the cell
- Fluxes computed using high-speed cell air temperature (Tcell) are compared to those computed using slow block temperature (Tblock)

SITES

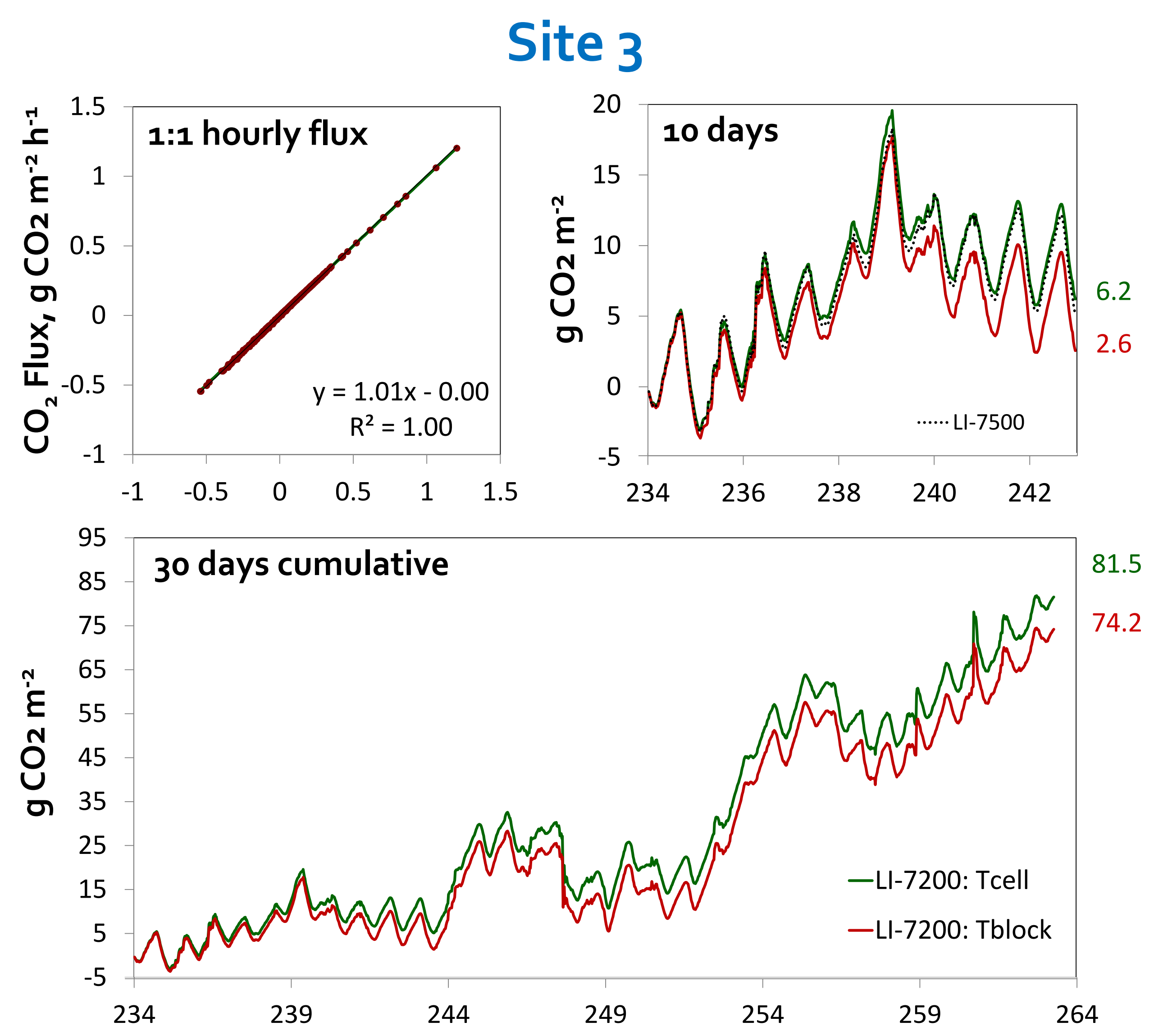
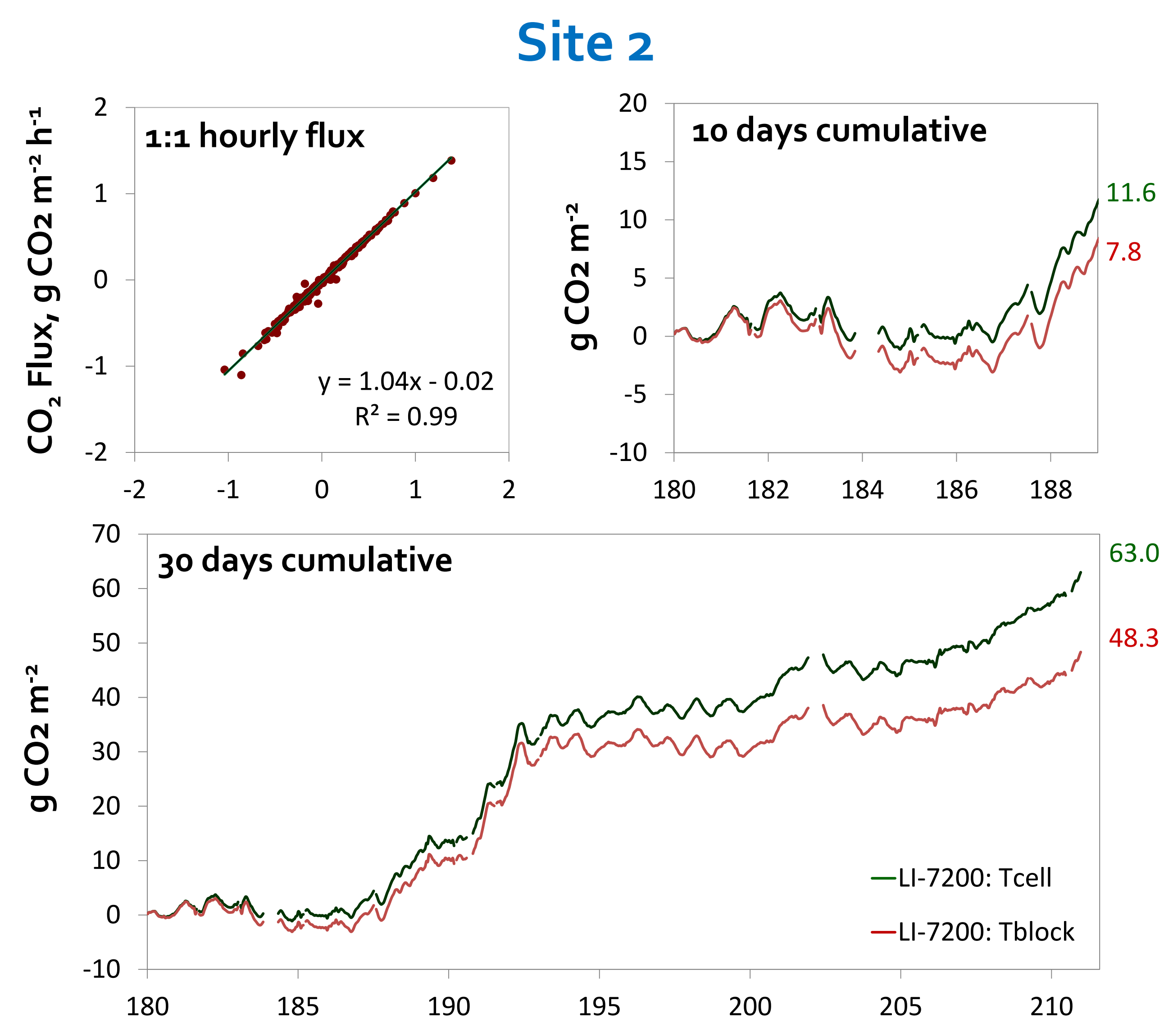
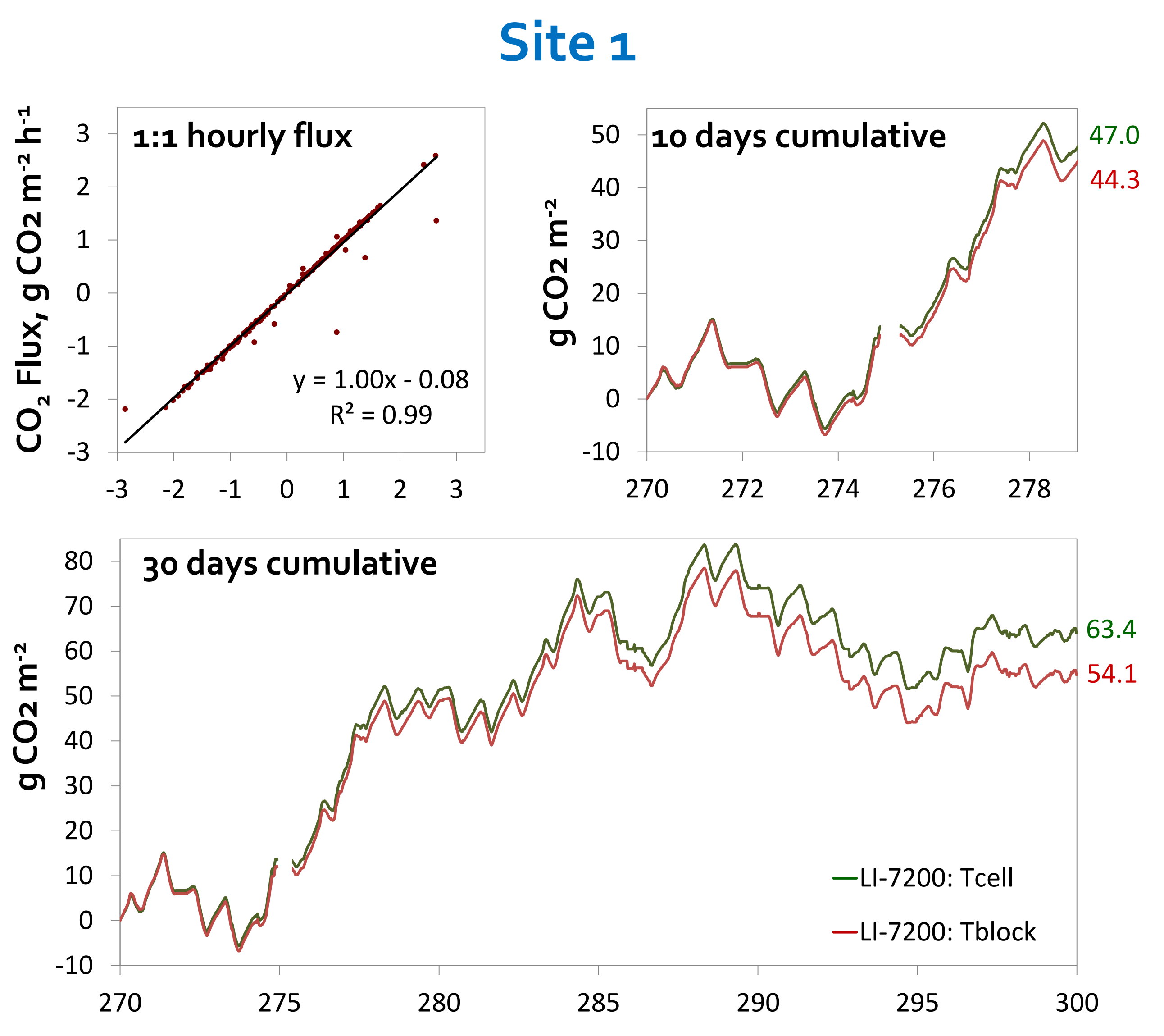
- Site 1:** LI-COR Experimental Research Station
Irrigated short grass, 0.05 m
3.8 m measurement height
Gill R350 + LI-7200 + EddyPro
- Site 2:** Himmelmoor site, U of Hamburg, Germany
Short grass, 0.05 m
8.0 m measurement height
Gill R3 + LI-7200 + EddyPro
- Site 3:** Temporary plot with ryegrass and weeds
Medium variable canopy, 0.1-0.6 m
2.5 m measurement height
CSAT3 + LI-7200 + EdiRe

SAMPLED AIR TEMPERATURE

- Example from Site 1
- Tcell and Tblock are different in means and fluctuations
- The differences are significant
- Spectra of Tcell and Tblock are shown
- Tube does not fully dampen fluctuations
- Residual fluctuations are captured by Tcell but not by Tblock



FLUX ERRORS RESULTED FROM USING SLOW TEMPERATURE



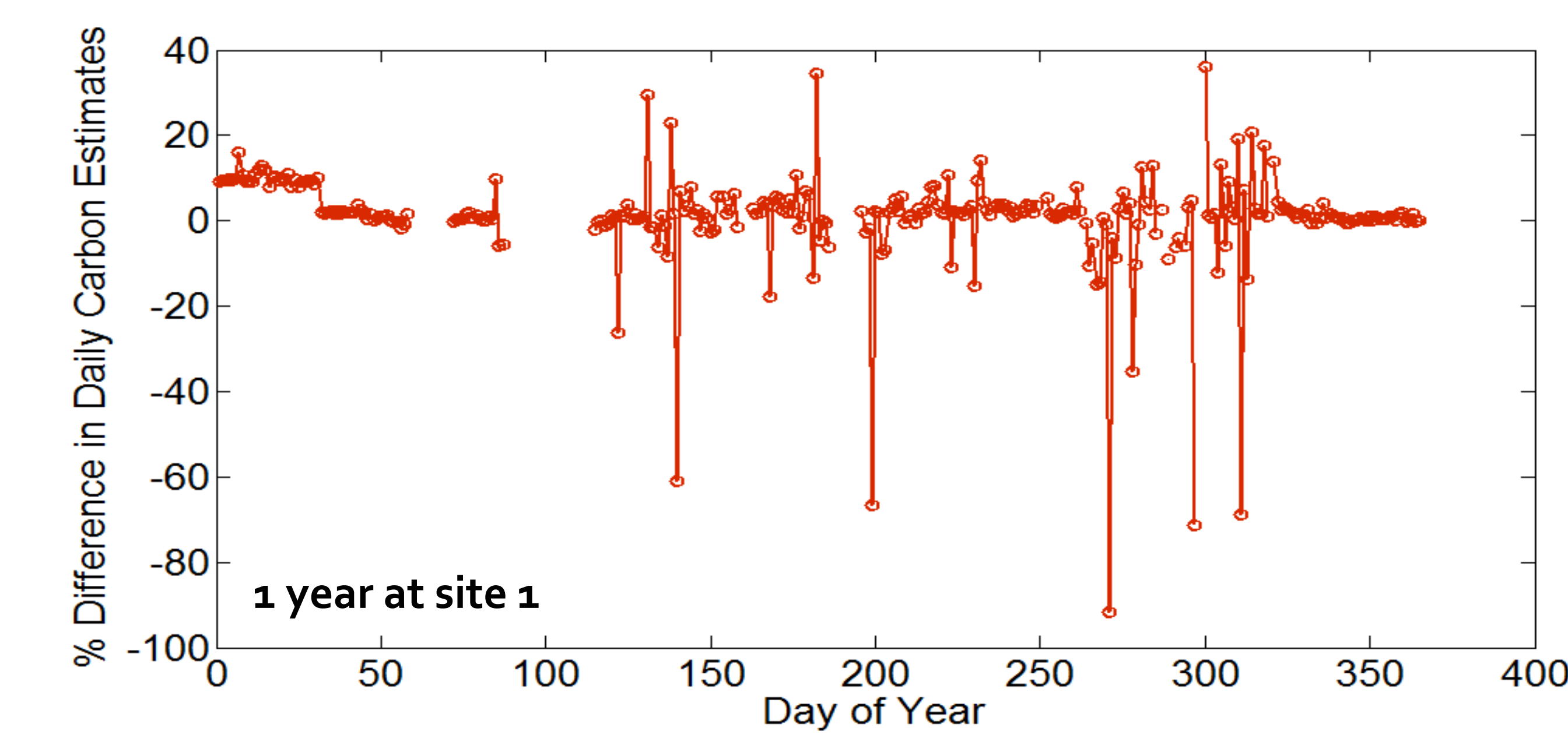
- The hourly 1:1 plot is deceptive, implying similar CO₂ fluxes when using fast Tcell and slow Tblock for calculating fast dry mole fraction.
- However, lack of fast air temperature measurements in the cell leads to accumulated error in CO₂ fluxes, noticeable after first 5 days and reaching 6% by the end of a 10-day period (47.0 vs 44.3 g CO₂ m⁻²). Over 30 fall days, the error increased to 9.3 g CO₂ (15% of monthly budget).

- Similar to site 1, the relatively close match of the 1:1 plot is deceptive (just 4% systematic flux underestimation from), but leads to a rapid accumulation of error in CO₂ flux, noticeable after first day, and significant by the end of the 10-day period (11.6 vs 7.8 g CO₂ m⁻²).
- The resulted error over 10 days is 3.8 g CO₂, or 33% of CO₂ budget. Over 30 summer days, the error increased to 14.7 g CO₂ (23% of monthly budget).

- At site 3, fluxes from LI-7200 computed using fast Tcell match open-path fluxes from LI-7500 computed using full WPL density terms
- The lack of fast air temperature leads to errors noticeable after first 3 days, and very significant by the end of the 10-day period (6.2 vs 2.6 g CO₂ m⁻²).
- The error over 10 days is 3.6 g CO₂, or 58% of the CO₂ budget. Over 30 days of senescence, the error increased to 7.3 g CO₂ (9% of monthly budget).

CONCLUSIONS

- Traditional analyzers used in eddy covariance have to use intake tubes 1000 times longer than the tube diameter to dampen most of ambient temperature fluctuation [2]
- In recently developed short-tube gas analyzers, fast temperature of sampled air must be measured inside the cell to avoid significant errors in CO₂ fluxes and budgets
- When cell block temperature is used instead of properly weighed fast cell air temperature, errors in CO₂ fluxes and carbon budgets can reach 60-90% on specific days, and can be as high as 58% over 10-day period
- Such error are expected to be significant in any environment but particularly detrimental for the sites where 24-hour or yearly budgets have comparable photosynthetic and ecosystem respiration components



REFERENCES

[1] Burba, G., D. McDermitt, D. Anderson, M. Furtaw, and R. Eckles RD. 2010. Novel design of an enclosed CO₂/H₂O gas analyser for eddy covariance flux measurements. *Tellus B*, 62, 743-748. [2] Rannik, U., Vesala, T., and Keskinen, R., 1997. On the damping of temperature fluctuations in a circular tube relevant to the eddy-covariance technique. *J. Geophys. Res.*, 102 (D11), 12,789-12,794. [3] Burba, G., A. Schmidt, R. Scott, T. Nakai, J. Kathilankal, G. Fratini, C. Hanson, B. Law, D. McDermitt, R. Eckles, M. Furtaw, and M. Velgersdyk, 2012. Calculating CO₂ and H₂O eddy covariance fluxes from an enclosed gas analyzer using an instantaneous mixing ratio. *Global Change Biology*, 18(1): 385-399

ACKNOWLEDGEMENTS

The flux measurements at the Himmelmoor site are run by the group of Prof. Lars Kutzbach at the Institute of Soil Science, University of Hamburg, and are funded by the Cluster of Excellence "CI-SAP" (EXC177), by the German Research Foundation (DFG)