

DIRECT CONTINUOUS MEASUREMENTS OF METHANE EMISSIONS FROM A LANDFILL

Method, Station and Latest Results

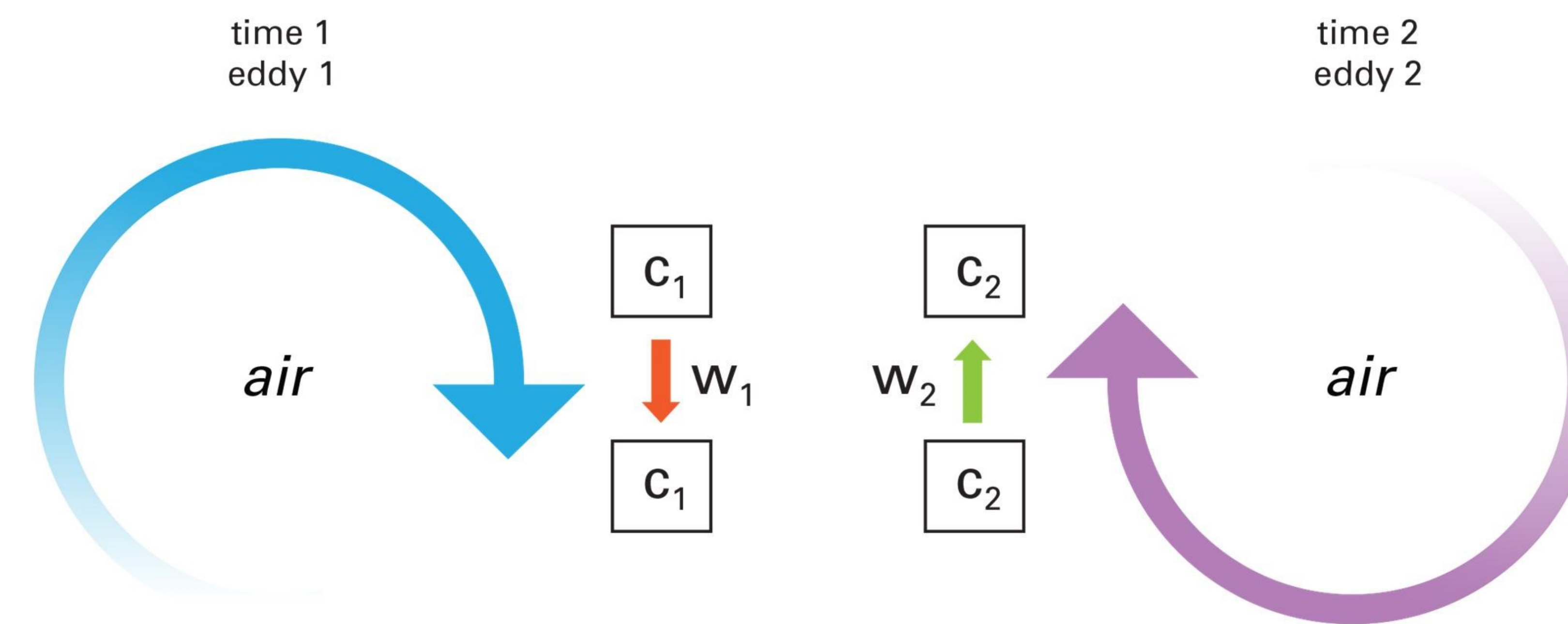
George Burba^{1*}, Liukang Xu¹, Xiaomao Lin², Jim Amen¹, Karla Welding³, and Dayle McDermitt¹

¹LI-COR Biosciences, Lincoln, NE, USA ²Kansas State University, Manhattan, KS, USA ³Bluff Road Landfill, Lincoln, NE, USA

INTRODUCTION

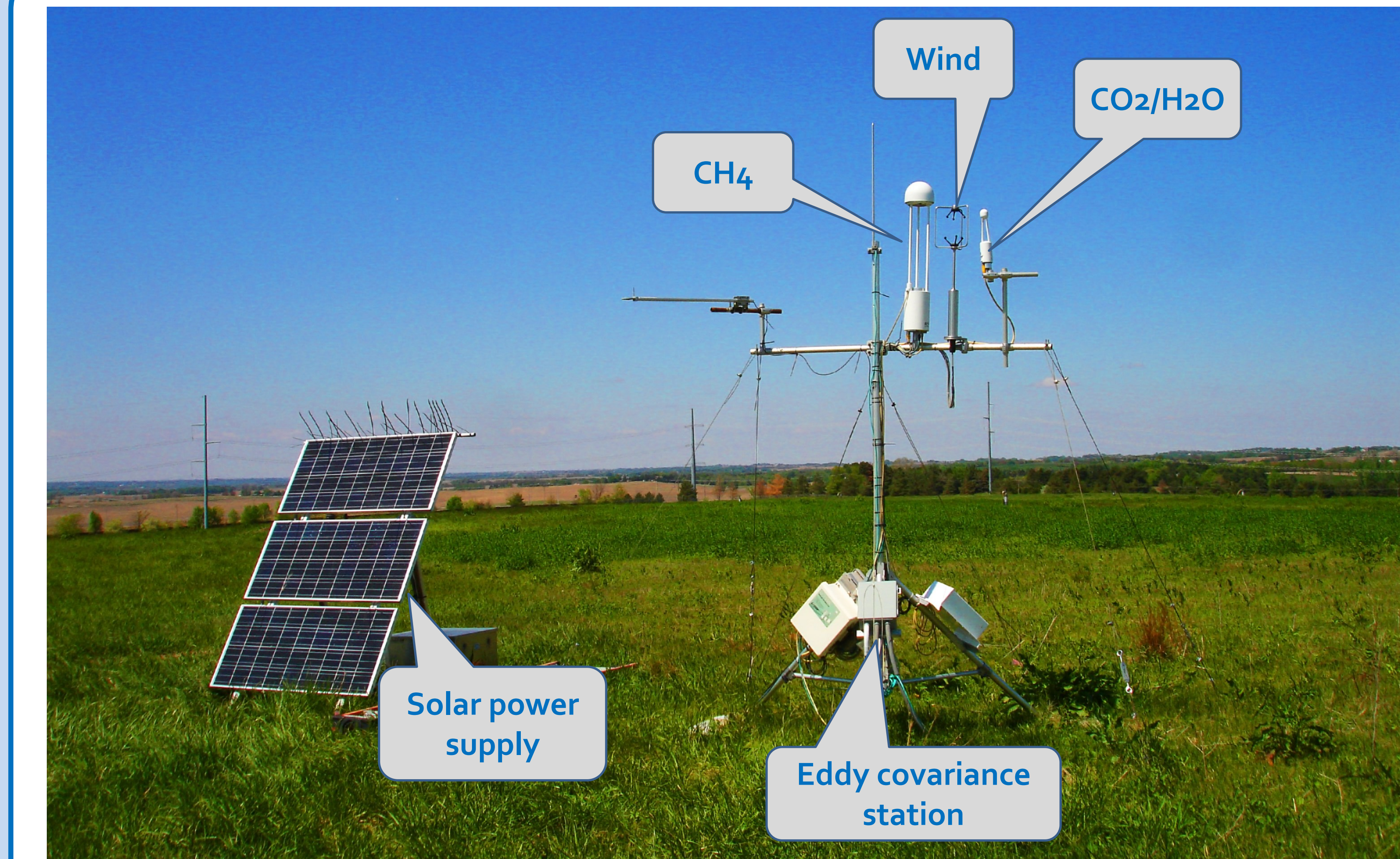
- Majority of year-round ecosystem-level CO₂, CH₄, and H₂O emissions measurements are presently done using eddy covariance method [1-3], with over 600 stations operating in over 120 countries, providing:
 - accurate measurements with high temporal resolution
 - time scales from hourly to daily and multi-year periods
 - detection of any rapid changes due to weather or human events
 - simultaneous detection of seasonal and yearly dynamics
- These features can be very helpful when investigating CH₄ emissions from landfills which historically have been greatly understudied [4-8]:
 - highly dynamic nature of actively filled landfill
 - intermittent and indirect nature of typical landfill measurements
 - widely variable results and conclusions
- Here we describe latest results from an eddy covariance CH₄/CO₂/H₂O station continuously operating for over 4 years in the center of a city landfill near Lincoln, Nebraska [9]

METHOD



- Wind moves air across the surface, creates turbulent vortices, or eddies
- 3D anemometer measures wind and gas analyzer measures gas content
- Covariance of wind and gas content describes the emission or flux rate
- Such measurements from landfills can help answer the key questions:
 - how large are direct emissions?
 - how variable are the emissions?
 - why do the emissions change?

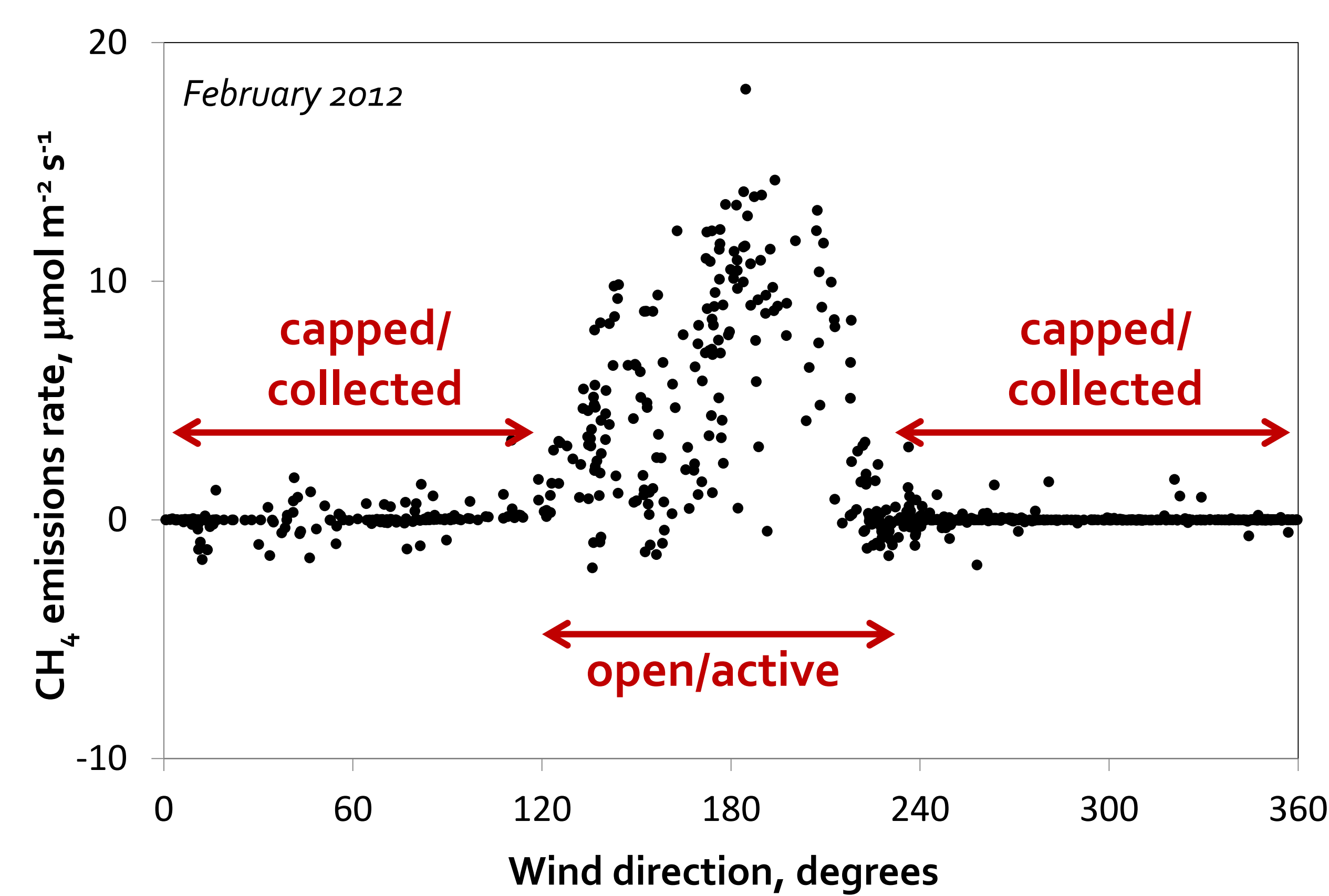
STATION



- Solar-powered wireless station positioned in the center of a landfill
- Automated online flux processing and quality control are used [10]
- Full access to raw data and final fluxes is provided from the office
- This approach helps gather novel results rapidly and with relative ease

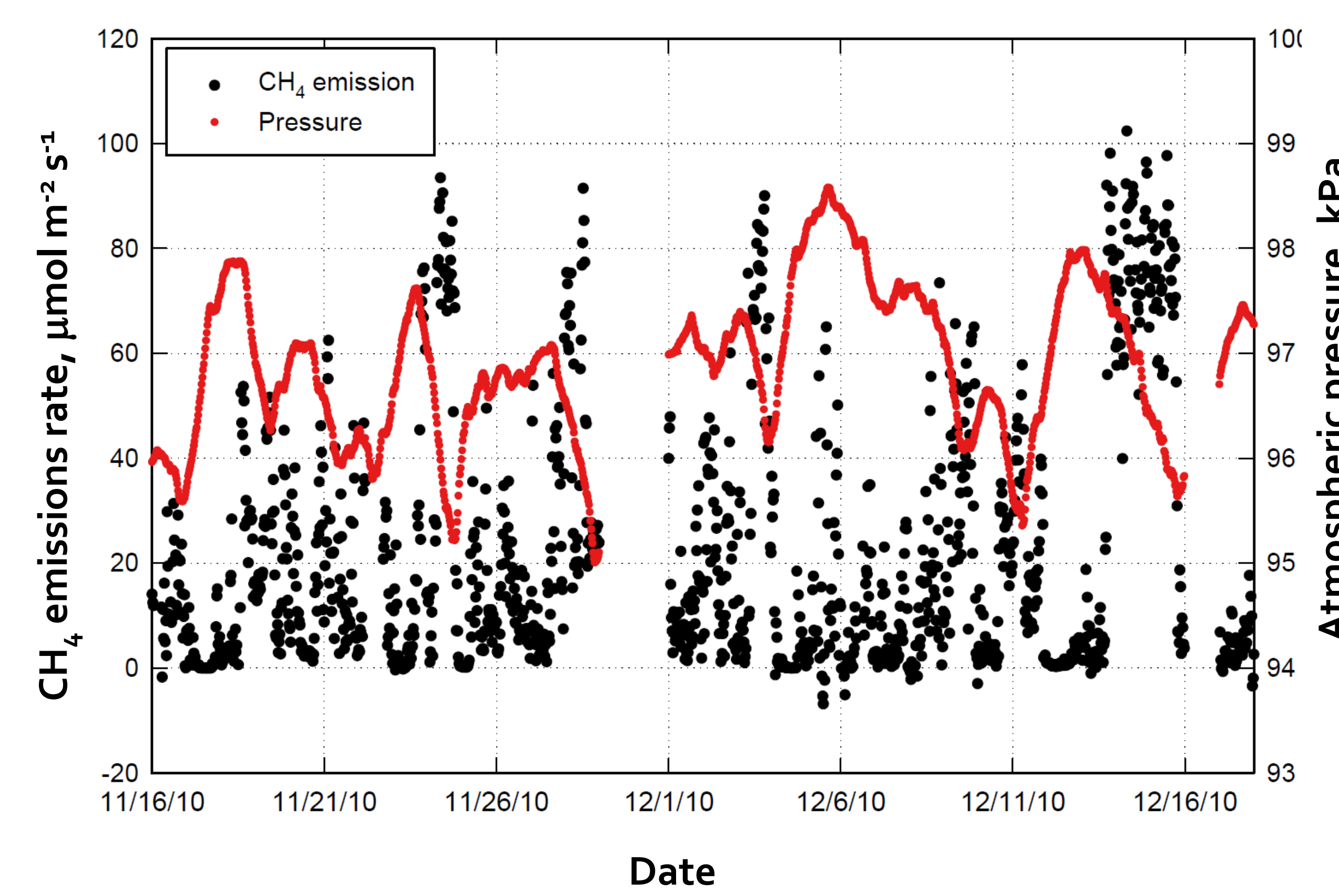
LATEST RESULTS

Footprint Coverage



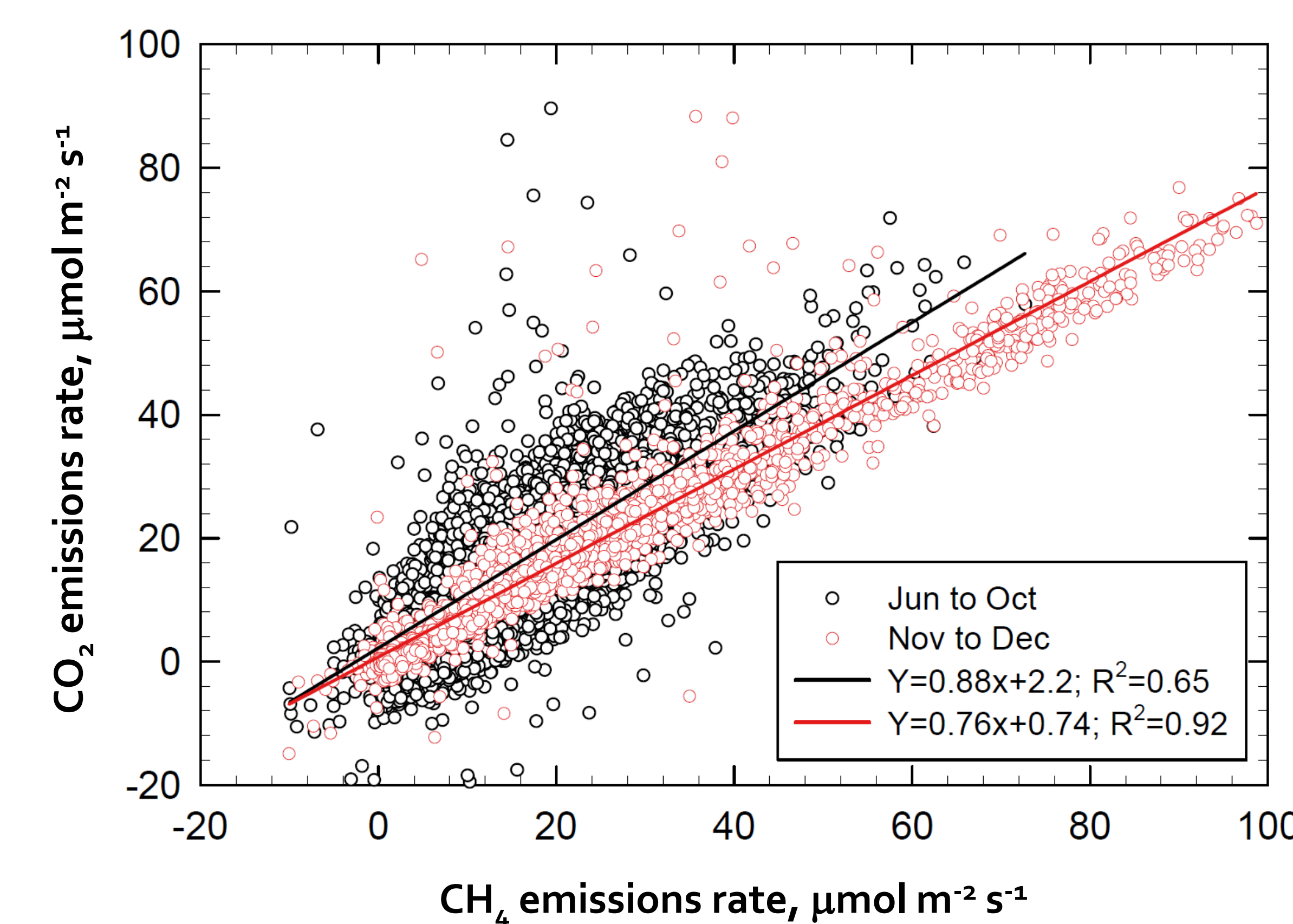
- Natural shifts in the wind direction allow effectively "scan" different parts of landfill at different times
- Effects of landfill capping are clearly evident when winds come from the capped area
- Fluxes from the capped area are reduced by orders of magnitude, from 10-20 µmol m⁻² s⁻¹ to near-zero

Pressure Effects



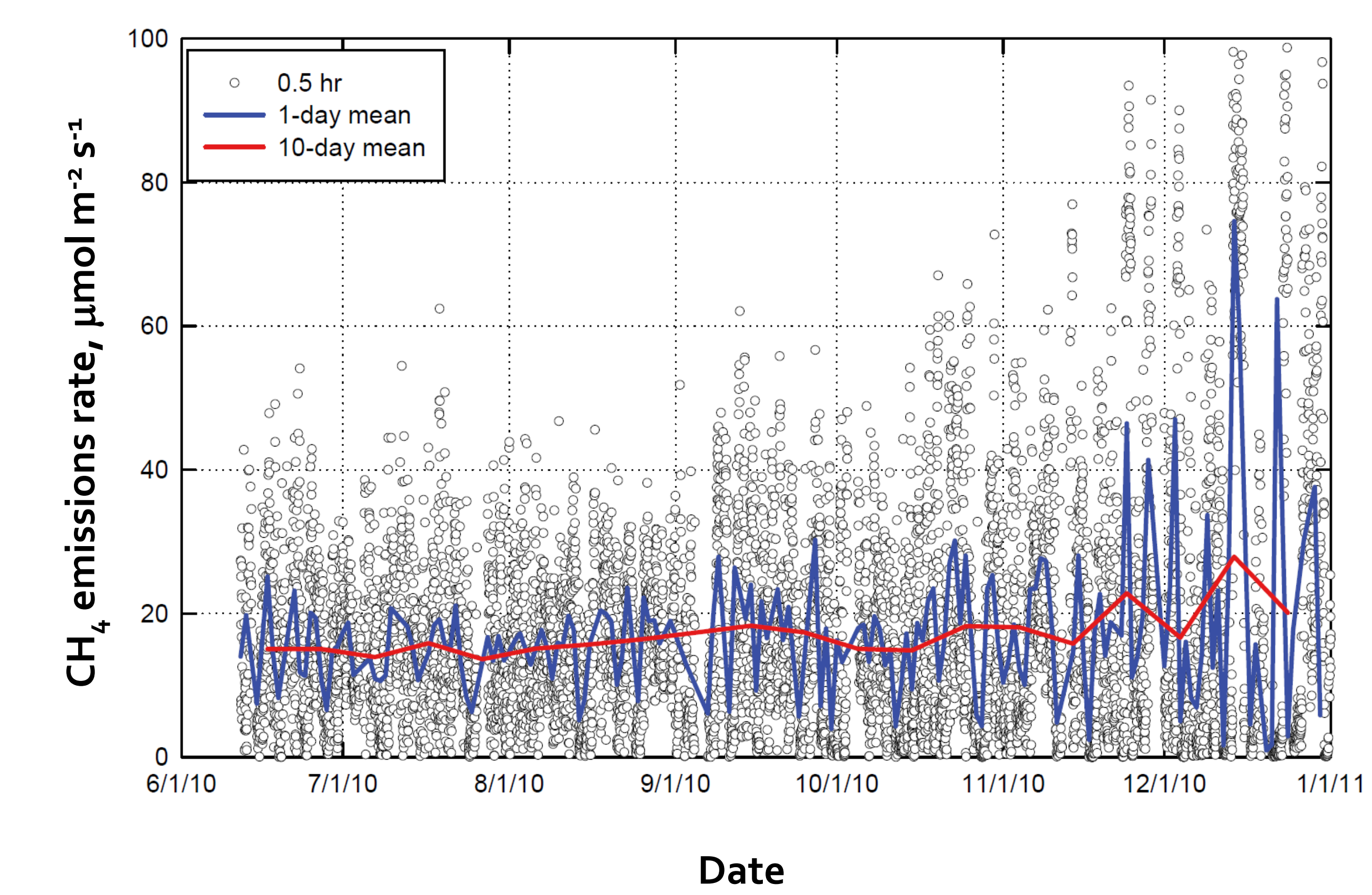
- Continuous measurements show what happens at variable atmospheric pressure
- Effects of changes in atmospheric pressure for this landfill is seen clearly in the plot above
- When pressure increases, fluxes are reduced orders of magnitude, from 60-100 µmol m⁻² s⁻¹ to near-zero

Oxidation Rates



- Concurrent CO₂ and CH₄ measurements allow a glimpse into oxidation process and rates
- Oxidation rates deduced from these data vary from near-zero in winter matching [11], to 7% in summer
- These results are specific to this landfill; results are quite different for a landfill in a different area [12]

Temporal Variability



- Continuous measurements show the extent of temporal variability in CH₄ landfill emissions
- At least 10 days of continuous measurements are needed for each state of weather or a season
- Occasional sporadic measurements of emitted CH₄ can easily be off by a few to tens orders of magnitude

REFERENCES

[1] Foken T. *Micrometeorology*. Springer-Verlag, 2008; [2] Aubinet M., Vesala T., Papale D. (Eds.). *Eddy Covariance: A Practical Guide to Measurement and Data Analysis*. Springer-Verlag, 2012; [3] Burba G. *Eddy Covariance Method for Scientific, Industrial, Agricultural and Regulatory Applications*. LI-COR Biosciences, 2013; [4] Bogner, J., M. Meadows, and P. Czepl. Fluxes of methane between landfills and the atmosphere: Natural and engineered controls. *Soil Use Management*, 1997 (13): 268-277; [5] Schroth, M. H., W. Eugster, K. E. Gómez, G. Gonzalez-Gil, P. A. Niklaus, and P. Oester. Above- and below-ground methane fluxes and methanotrophic activity in a landfill-cover soil. *Waste Management*, 2013 (32): 879-889; [6] Bingemer, H. G., and P. J. Crutzen. The production of methane from solid wastes. *J. Geophys. Res.*, 1987 (92): 2181-2187; [7] Czepl, P. W., B. Mosher, R. C. Harriss, J. H. Shorter, J. B. McManus, and C. E. Kolb. Landfill methane emissions measured by enclosure and atmospheric tracer methods. *J. Geophys. Res.*, 1996 (101): 16711-16719; [8] Giani, L., J. Bredenkamp, and I. Eden. Temporal and spatial variability of the CH₄ dynamics of landfill cover soils. *J. Plant Nutr. Soil Sci.*, 2002 (165): 205-210; [9] Xu, L., X. Lin, J. Amen, K. Welding, and D. McDermitt. Impact of changes in barometric pressure on landfill methane emission. *Global Biogeochem. Cycles*, 2014 (28): 17 pp.; [10] LI-COR Biosciences. Practical Solutions for Large-scale CO₂, H₂O, and CH₄ Emission Measurements. *Environmental Technology*, 2013 (2): 8-9; [11] Spokas, K., and J. Bogner. Limits and dynamics of methane oxidation in landfill cover soils. *Waste Management*, 2011 31(5):823-832; [12] LI-COR Group, et al. Quantifying total landfill methane emission: methodology comparison between eddy-covariance and plume-trace methods. *AWMA 108th Annual Conference*, 2015 (in progress).