

Unlocking the (Reversal) Potential of SSM Electrophysiology: Transporter Stoichiometry with the SURFE²R N1

Nathan Thomas

Henzler-Wildman Lab

University of Wisconsin – Madison



WISCONSIN
UNIVERSITY OF WISCONSIN-MADISON



Integrated Program in
BIOCHEMISTRY

Stoichiometry is a crucial determinant of transporter function

LacY

1 proton per lactose

(West and Mitchell, 1972)

PepT_{sa}

5 protons per dipeptide

(Parker, Mindell, and Newstead, 2014)

-160mV PMF can drive:

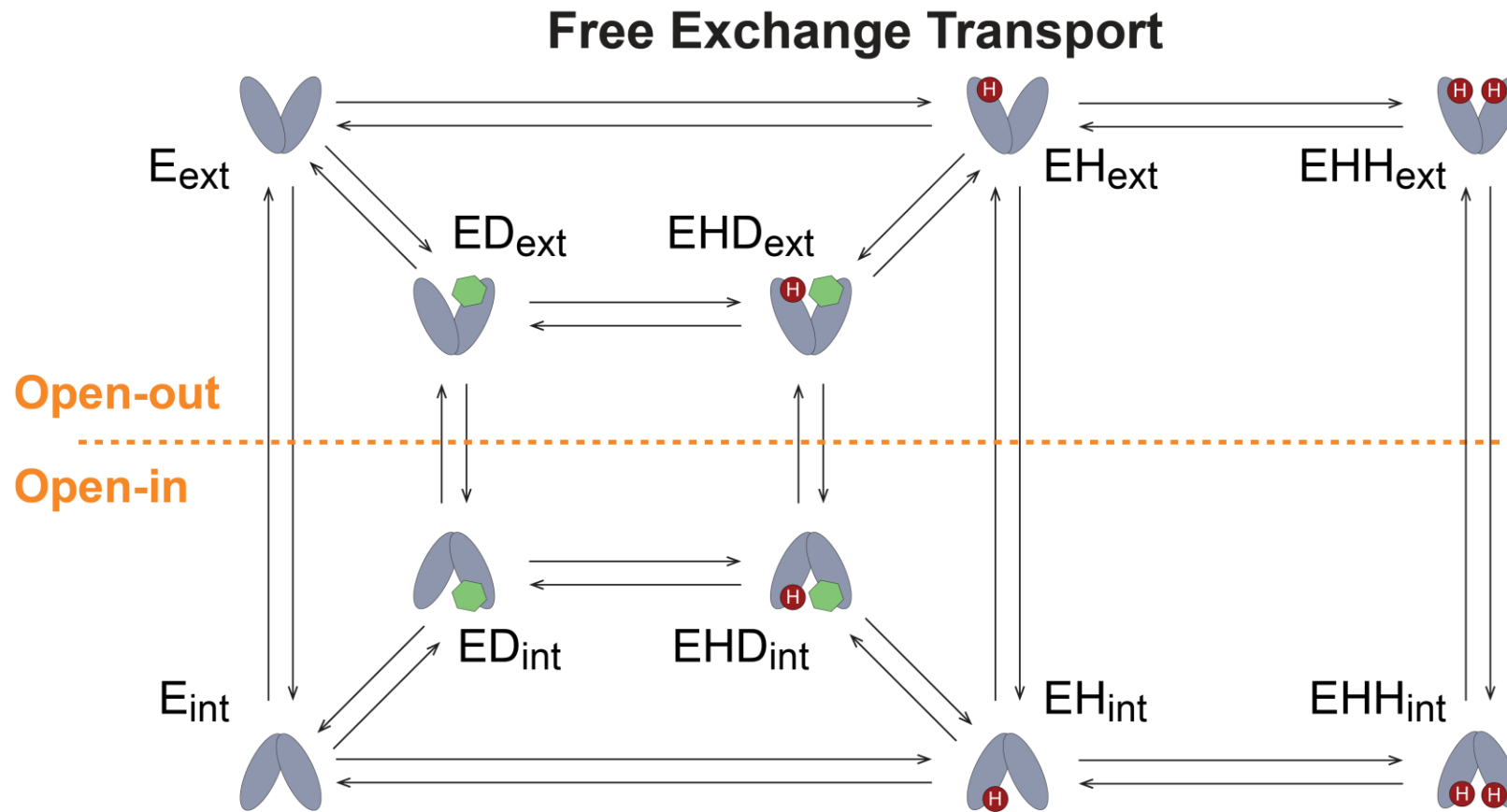
400-fold lactose gradient

-160mV PMF can drive:

10 trillion-fold dipeptide gradient

*Assumes tight coupling

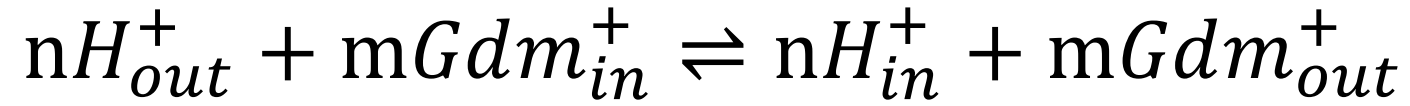
MDR efflux pump EmrE is not tightly coupled



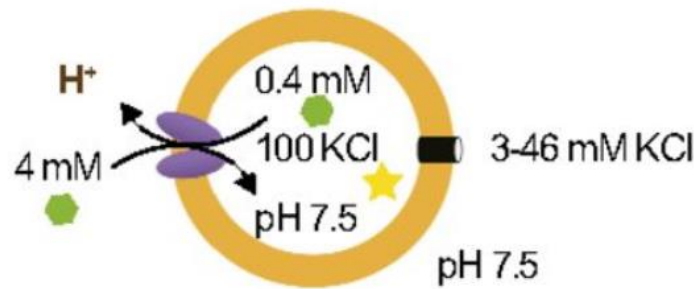
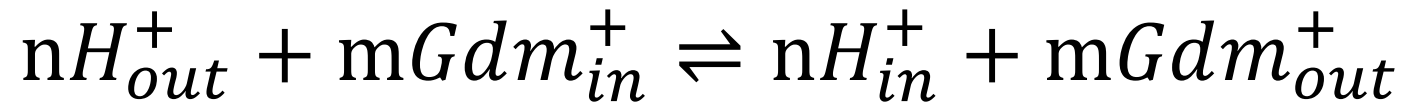
Multiple pathways – multiple stoichiometries?

How to measure transport stoichiometry

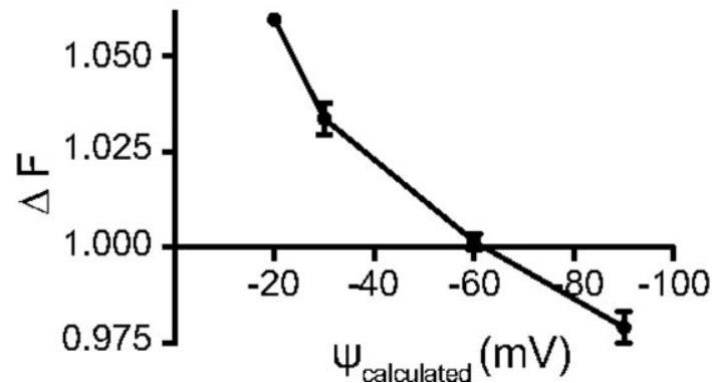
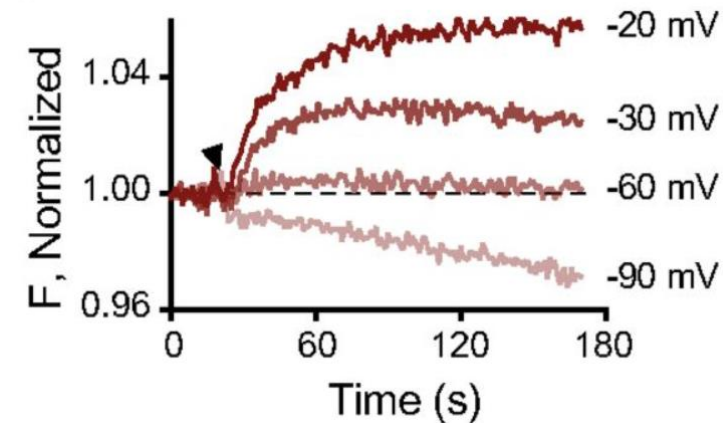
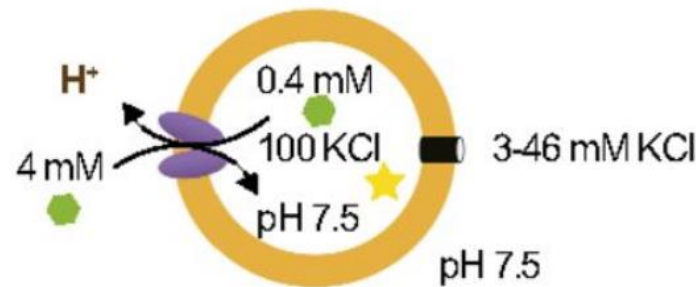
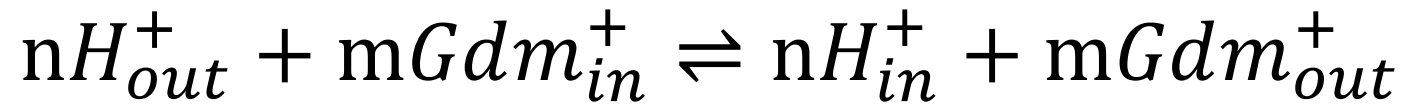
How to measure transport stoichiometry: Traditional Reversal Potential assay with GdX



How to measure transport stoichiometry: Traditional Reversal Potential assay with GdX

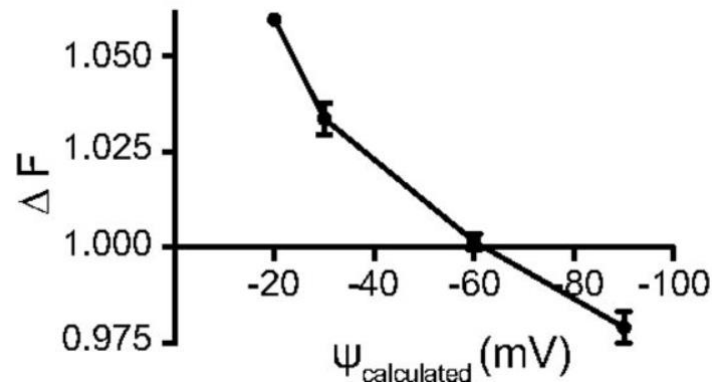
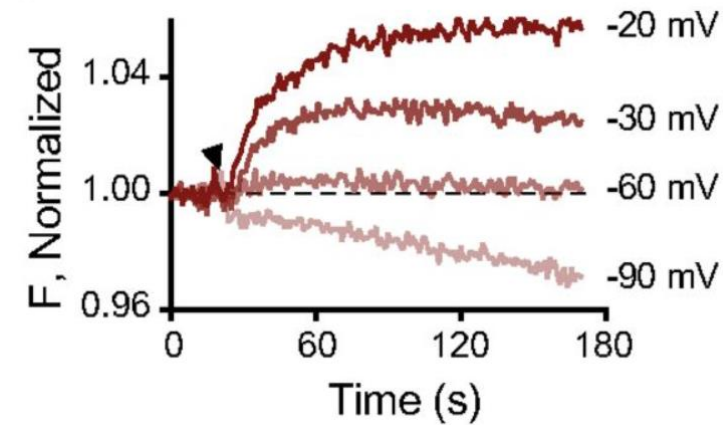
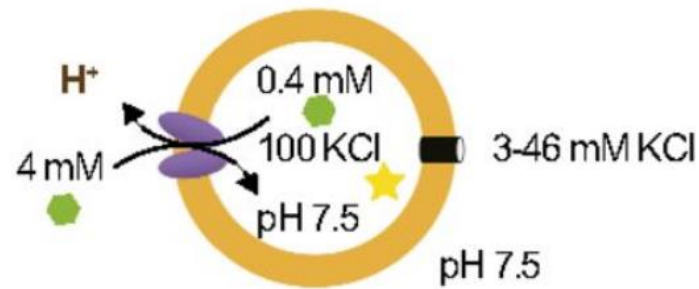
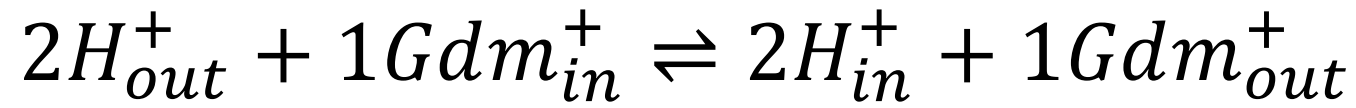


How to measure transport stoichiometry: Traditional Reversal Potential assay with GdX



$$E_{rev} = \frac{n\mu_i + m\mu_s}{-F(nz_i + mz_s)}$$

How to measure transport stoichiometry: Traditional Reversal Potential assay with GdX



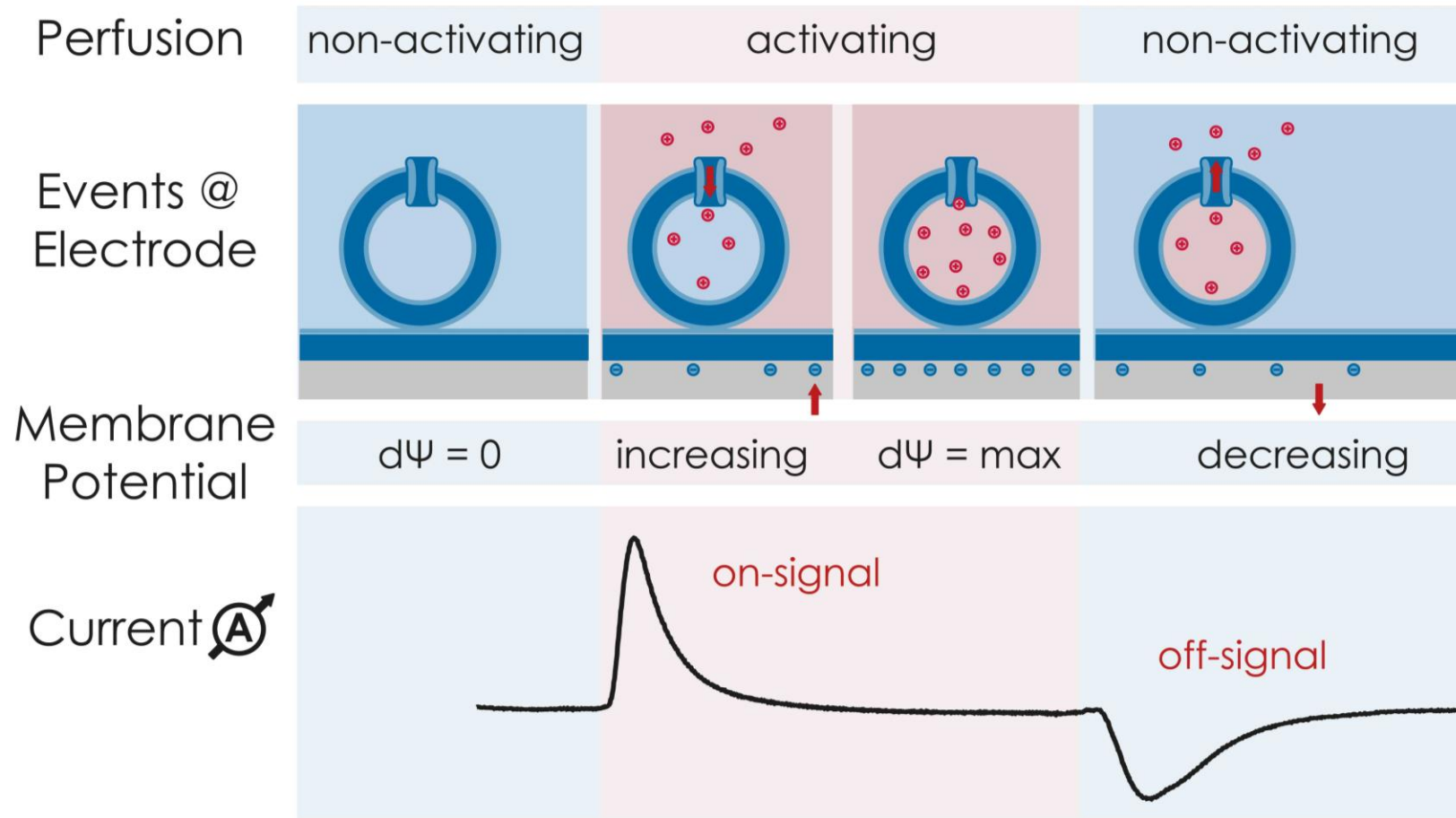
$$E_{rev} = \frac{n\mu_i + m\mu_s}{-F(nz_i + mz_s)}$$

Issues with reversal potential assays

- Technically difficult and time consuming
 - Need to ensure internal solutions are accurate
 - Different internal conditions require separate reconstitutions
- Need fluorophores or radioactivity to monitor transport
- Relatively large amounts of sample required

Can SSM-electrophysiology be
used to measure stoichiometry?

Overview of an SSM-electrophysiology experiment



SURFE²R N1

Reversal Potential Assay

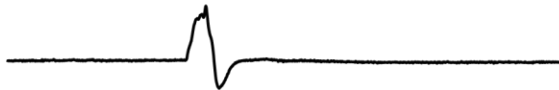
Assay Scheme – Proton gradient drives guanidinium transport

Buffer Flow:

100nM H⁺ —————→
1mM Gdm⁺



Current:



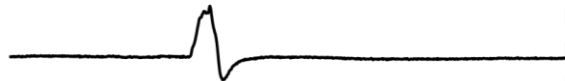
Assay Scheme – Guanidinium gradient drives proton transport

Buffer Flow:

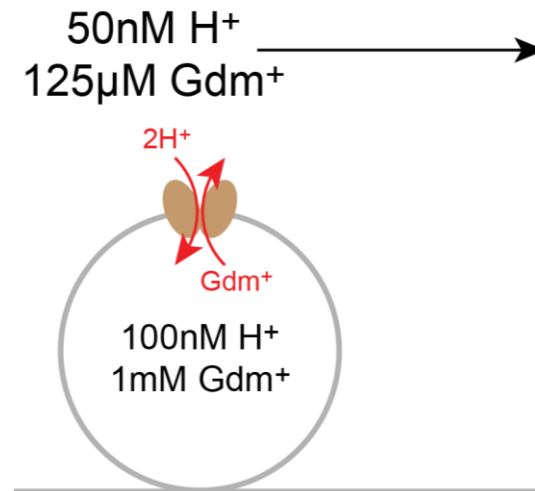
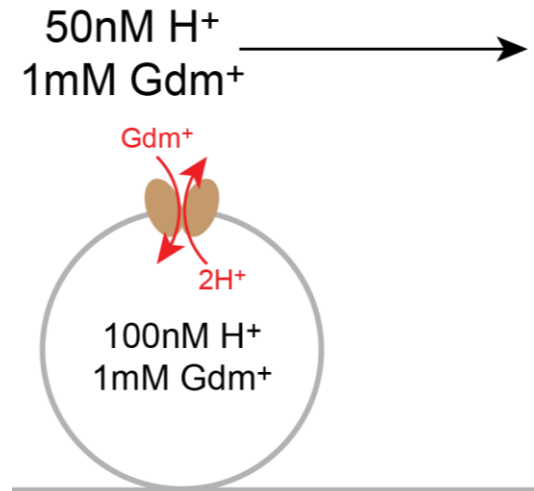
100nM H⁺
1mM Gdm⁺ →



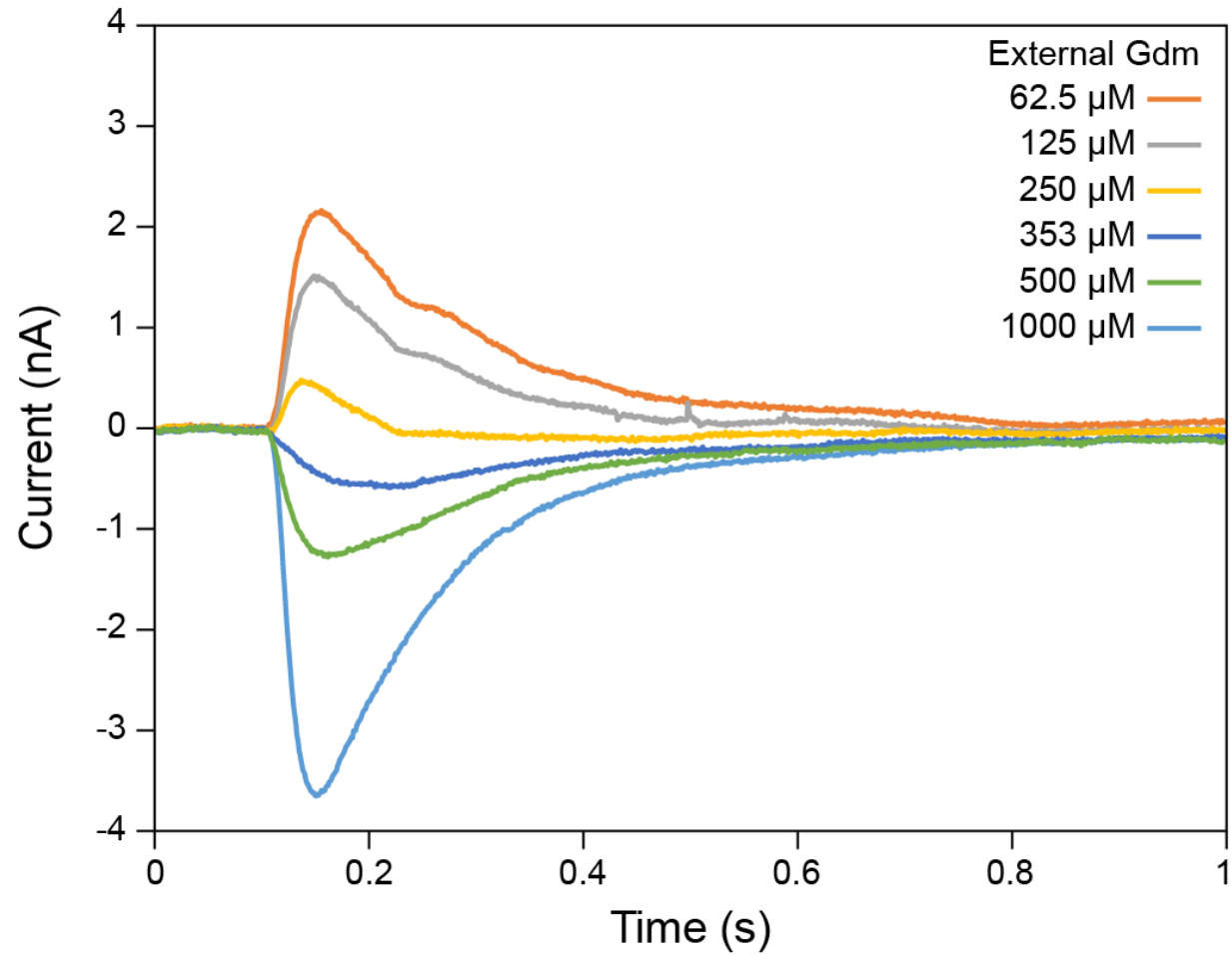
Current:



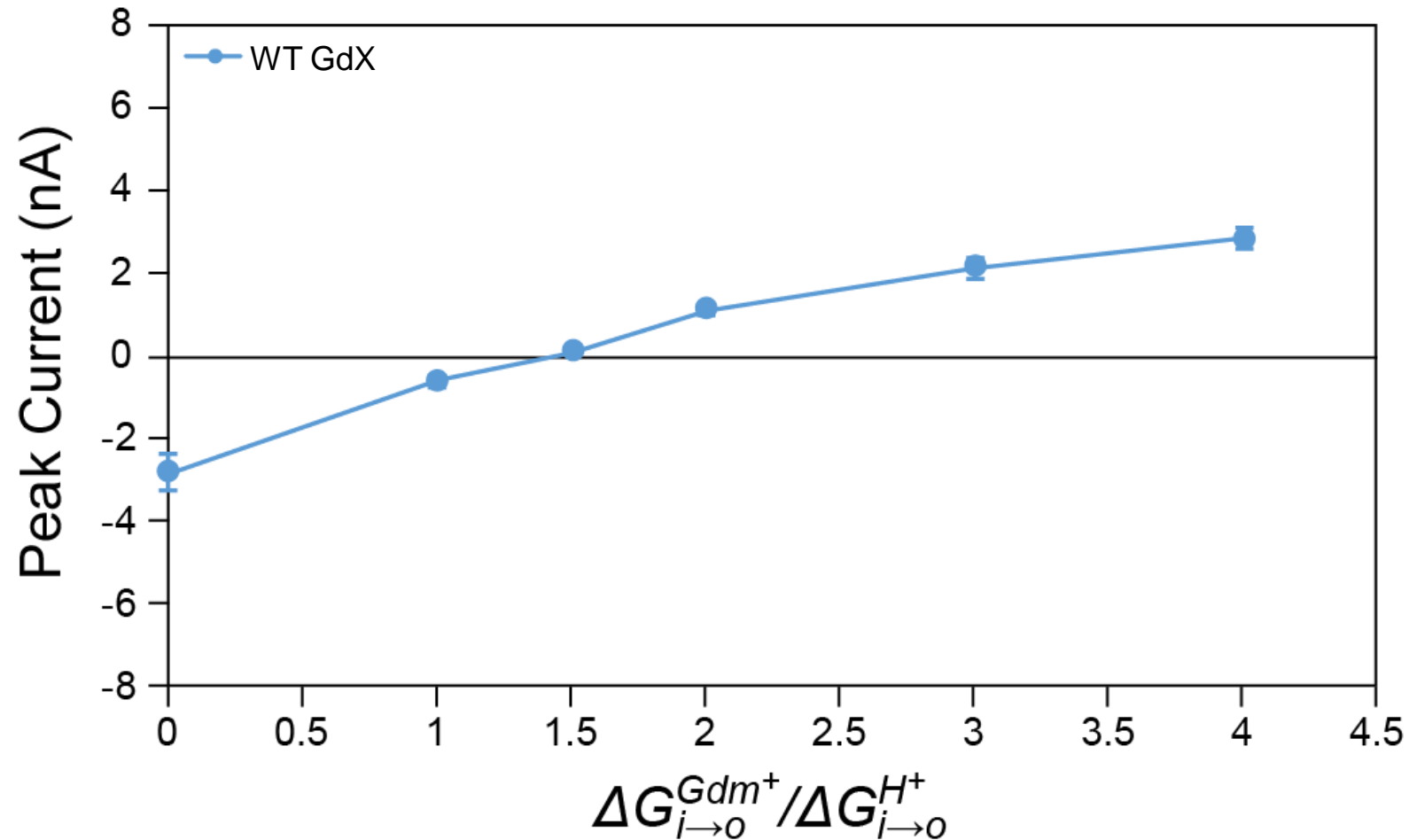
Transport current can be reversed



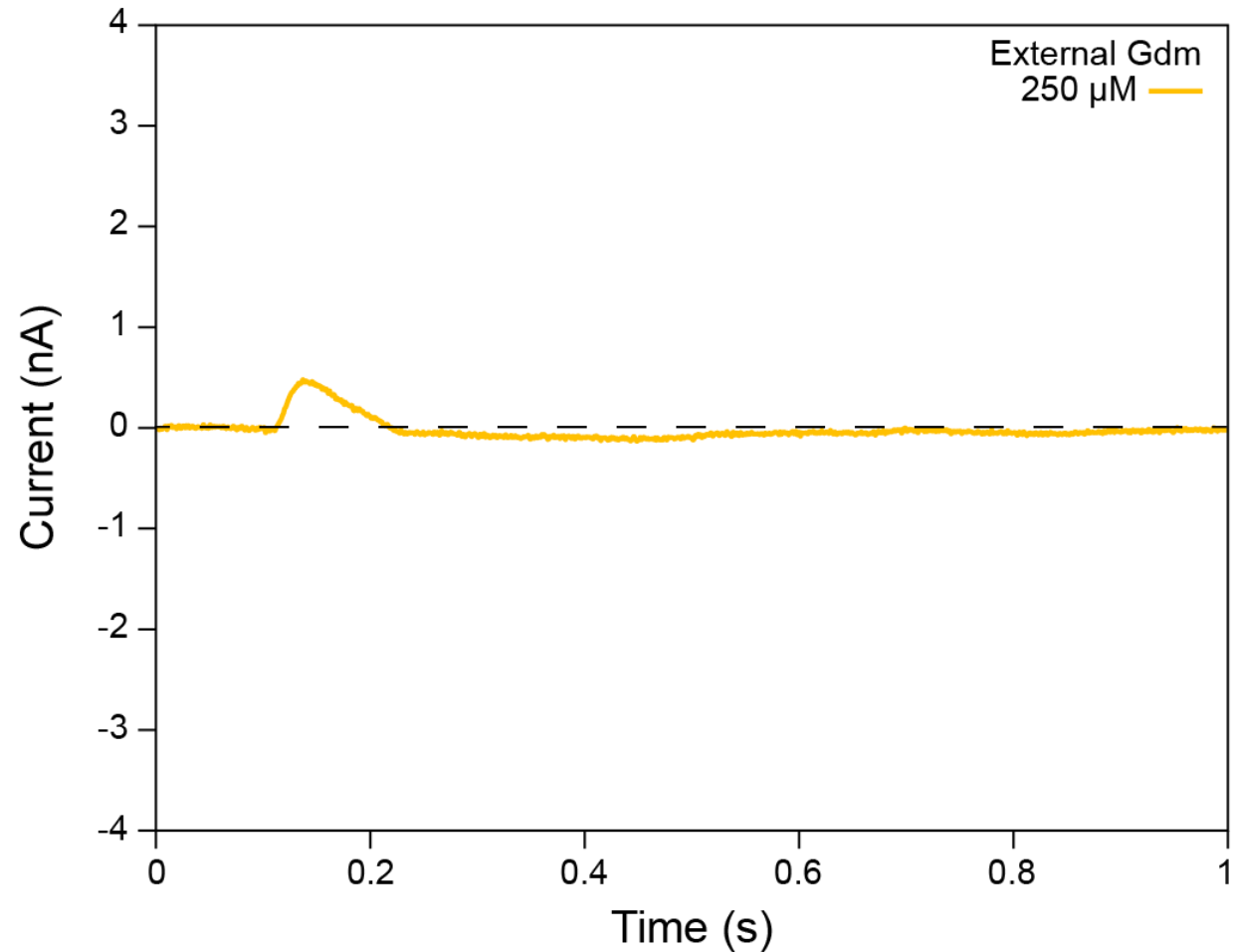
Transport current can be reversed



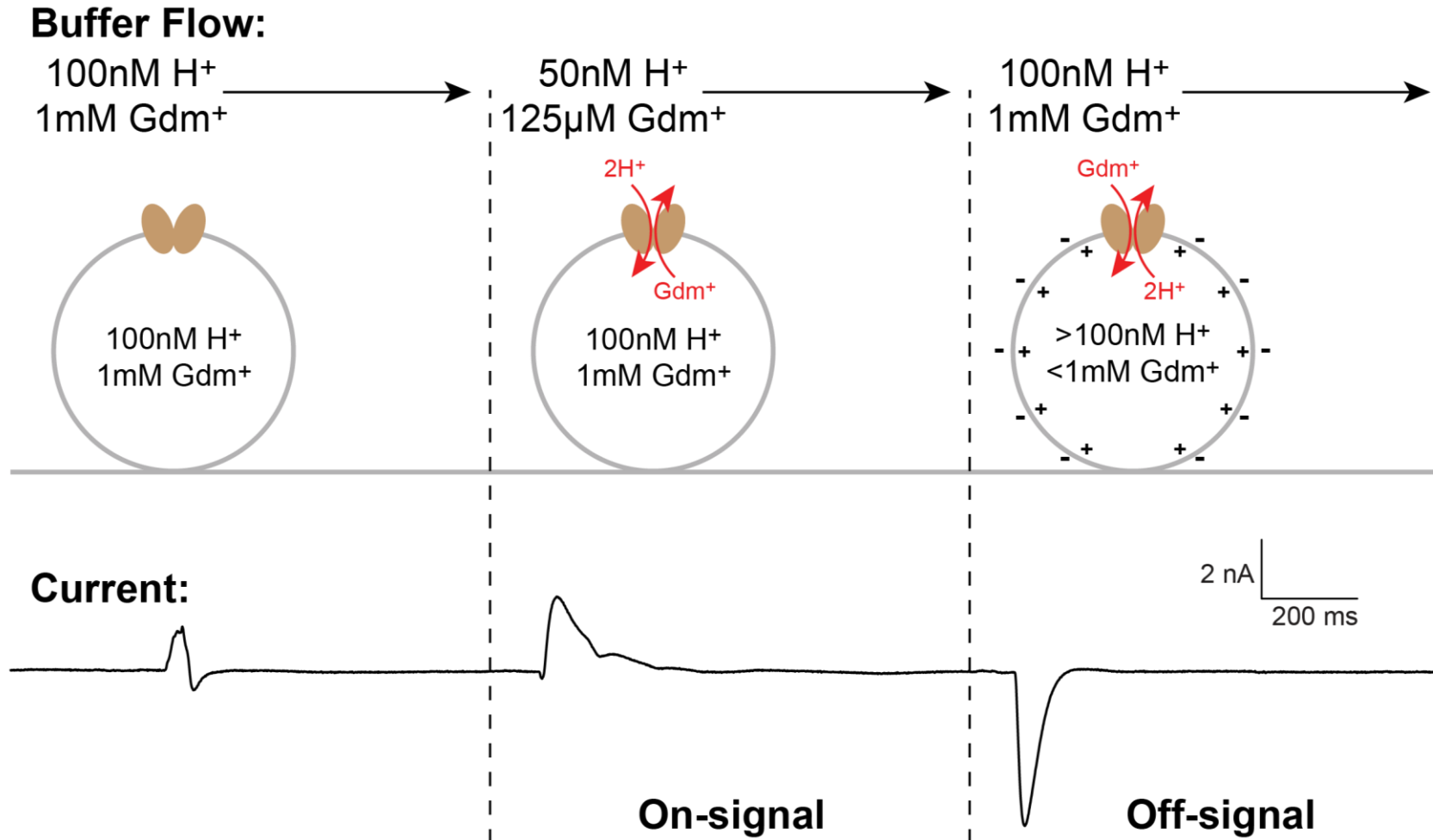
Peak current does not reverse at published reversal potential



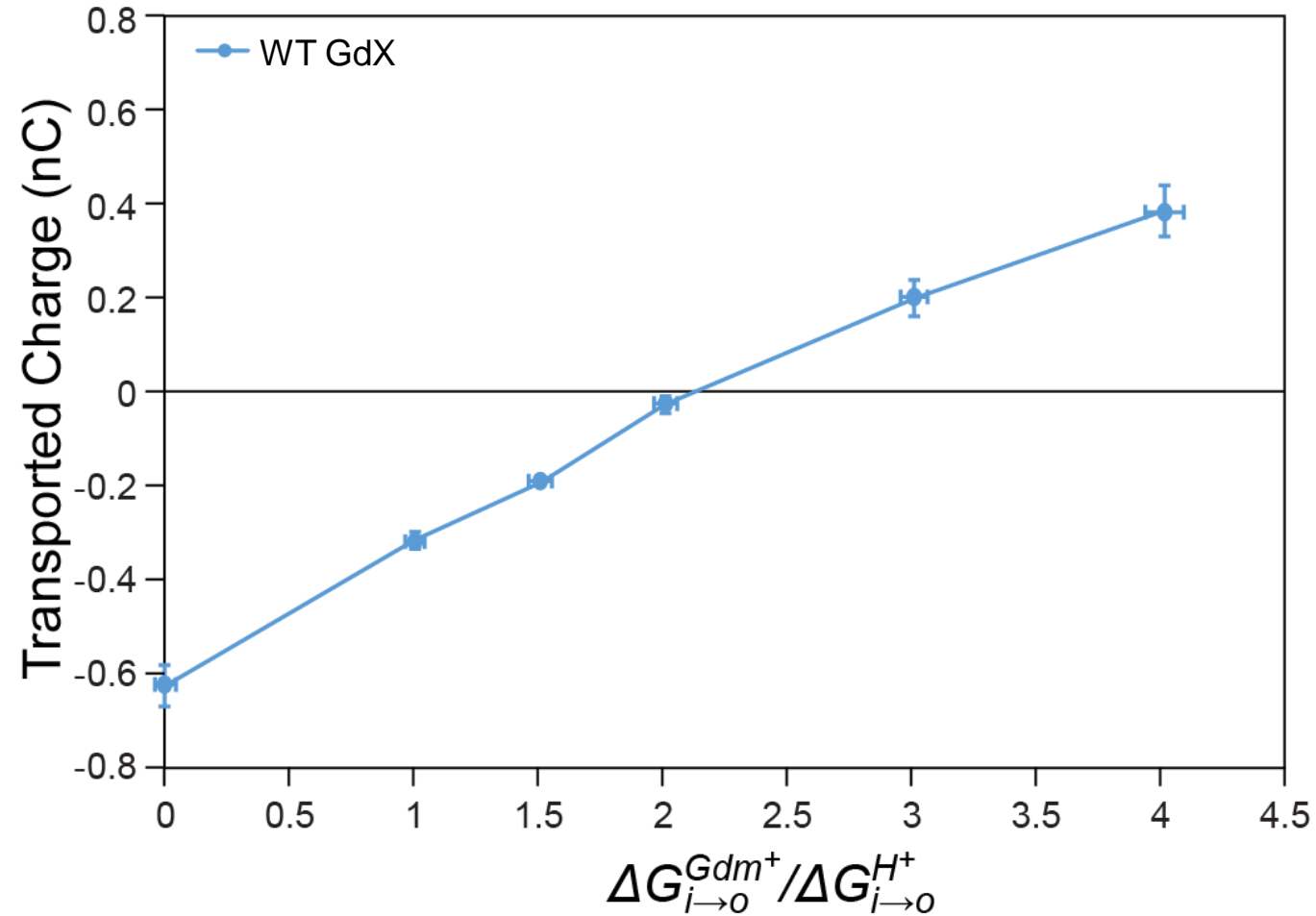
Peak current does not reverse at published reversal potential



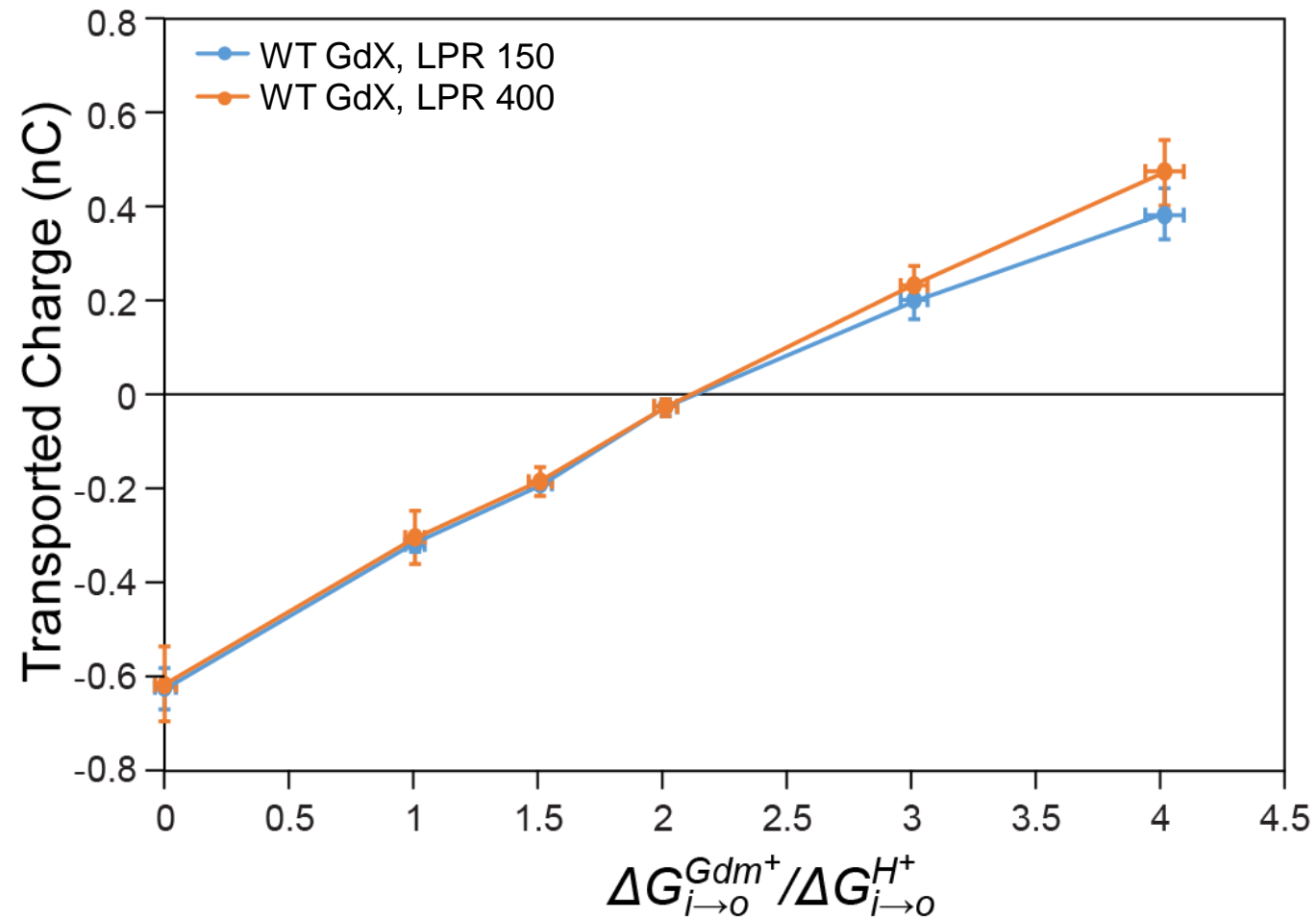
Integrate on-signal



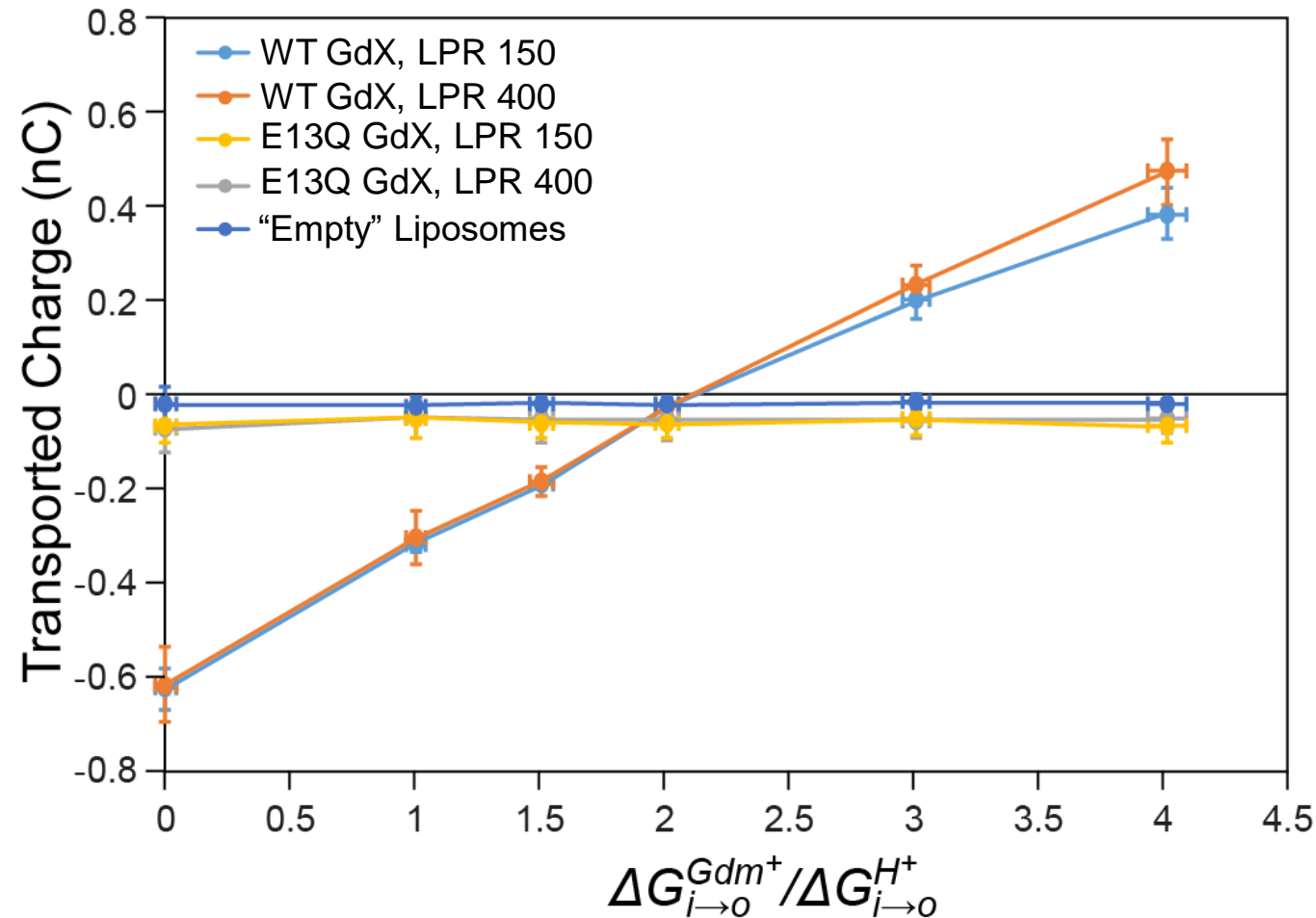
Integrated signal agrees with published stoichiometry



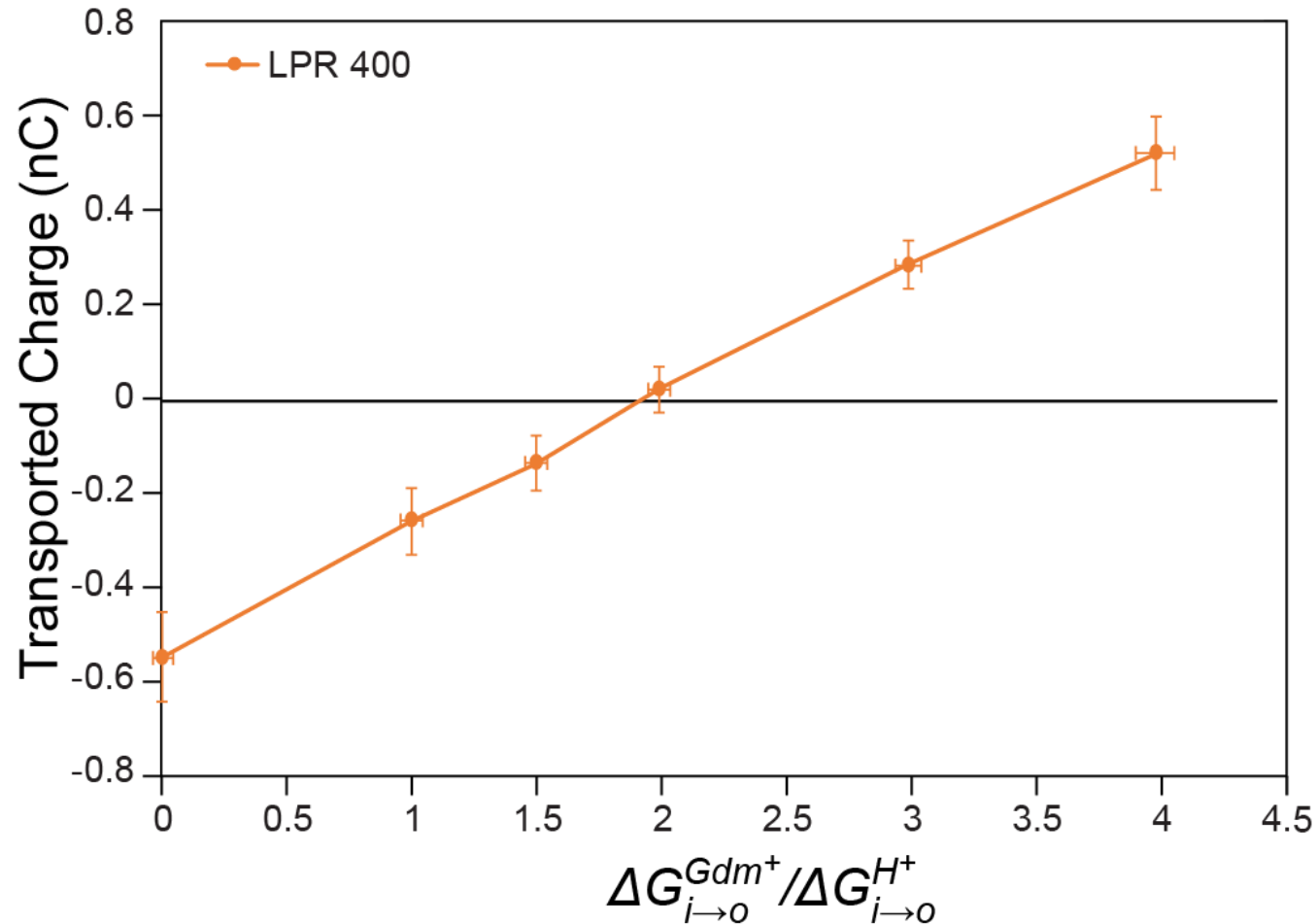
Signal is independent of protein concentration



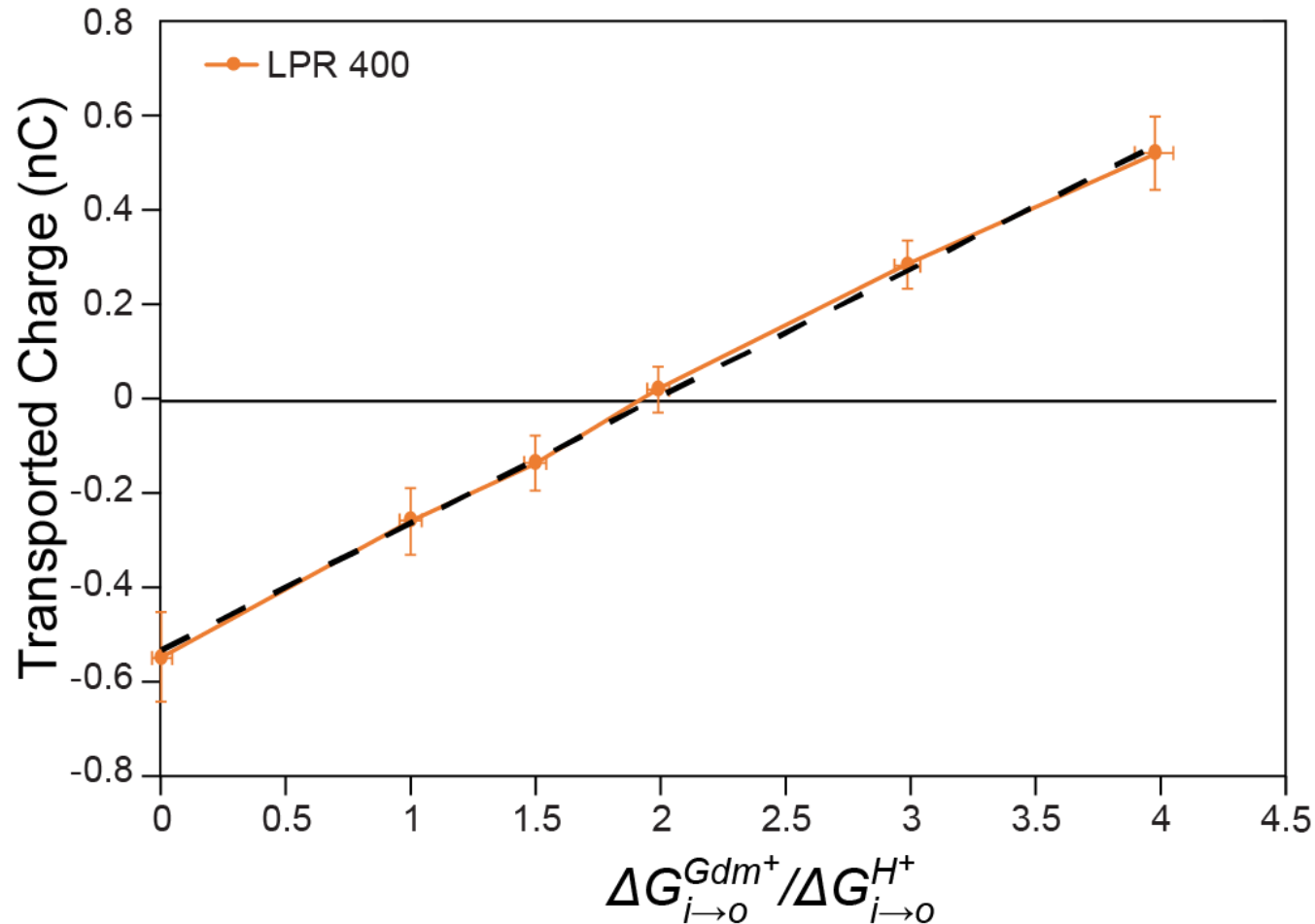
No significant currents for empty liposomes or transport-dead mutants



Transported Charge is a linear function of chemical potential



Transported Charge is a linear function of chemical potential



$$R^2 = 0.998$$

$$\text{x-intercept} = 1.98 \pm 0.06$$

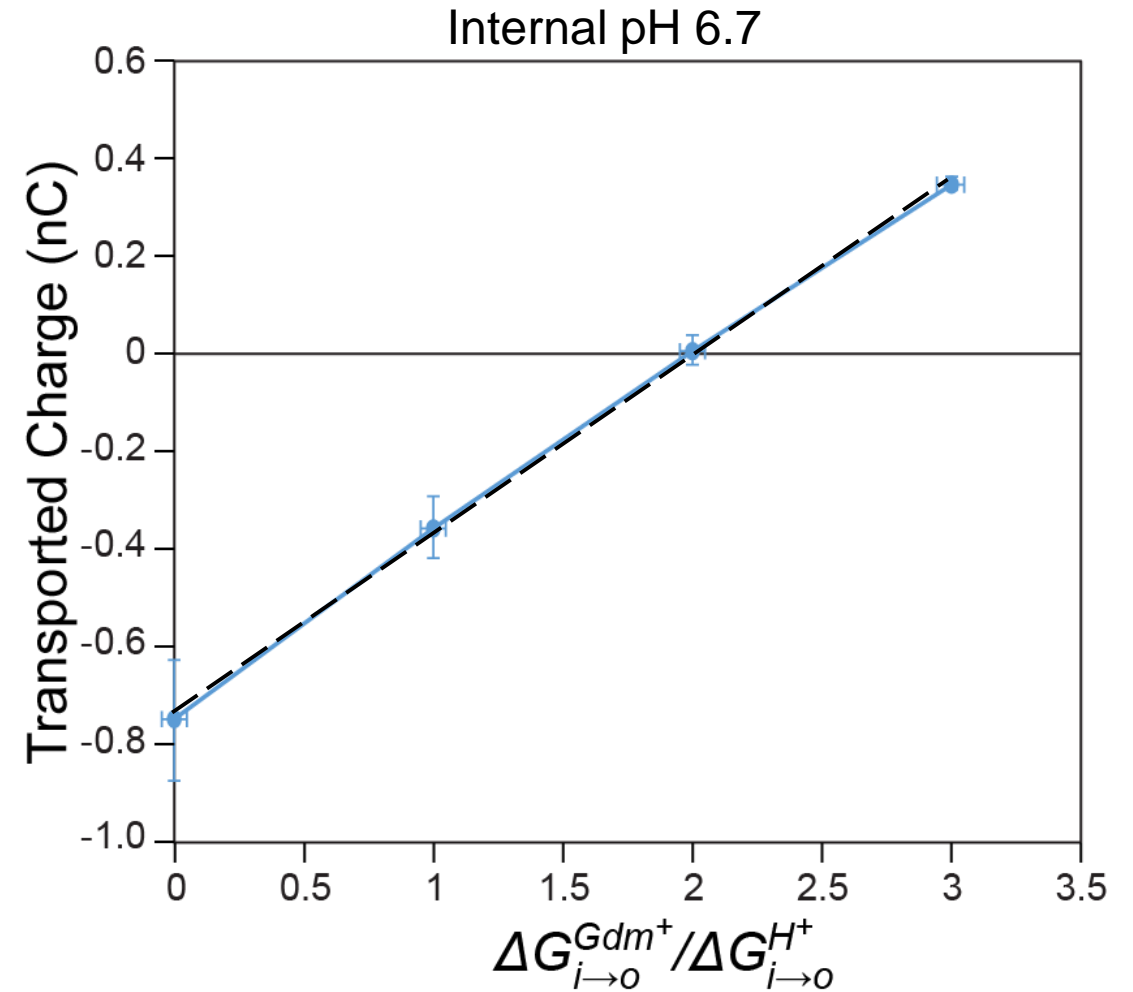
Sensor internal solution can be changed

- Reconstitute and prepare sensors at pH 7, 1 mM Gdm⁺
- Flow pH 6.7, 1 mM Gdm⁺ buffer over sensor
 - Monitor solution exchange until no current is observed – about 3 mL
- Repeat reversal potential assay
 - Internal buffer: 200 nM H⁺, 1 mM Gdm⁺
 - External buffer: 100 nM H⁺, varied Gdm⁺

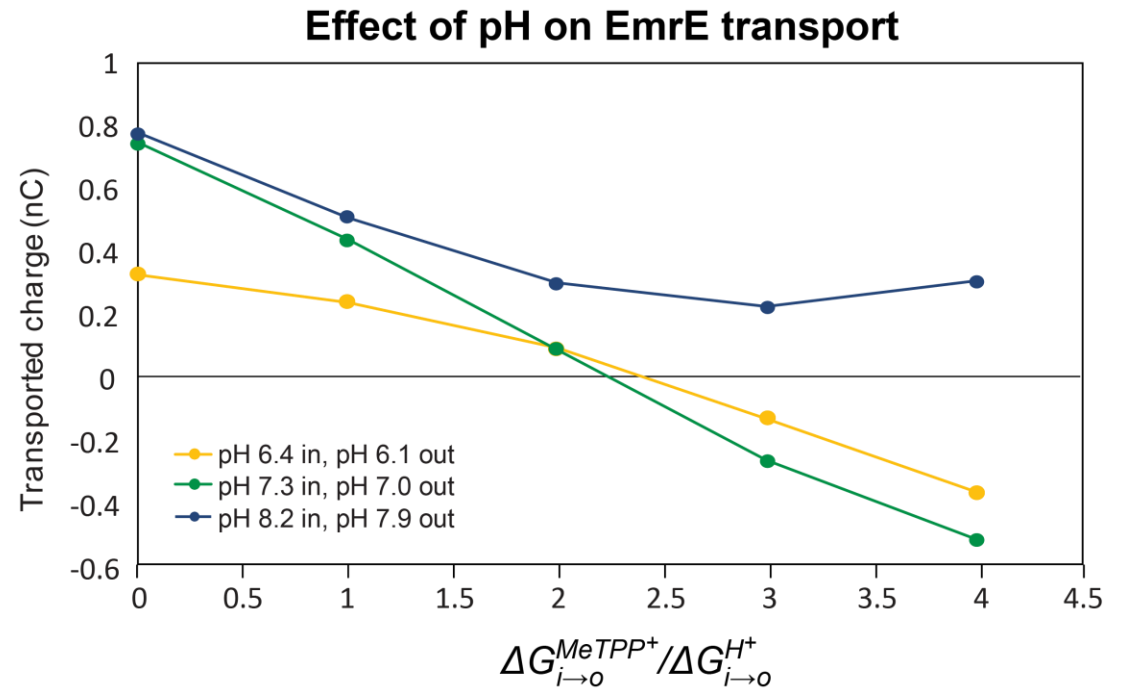
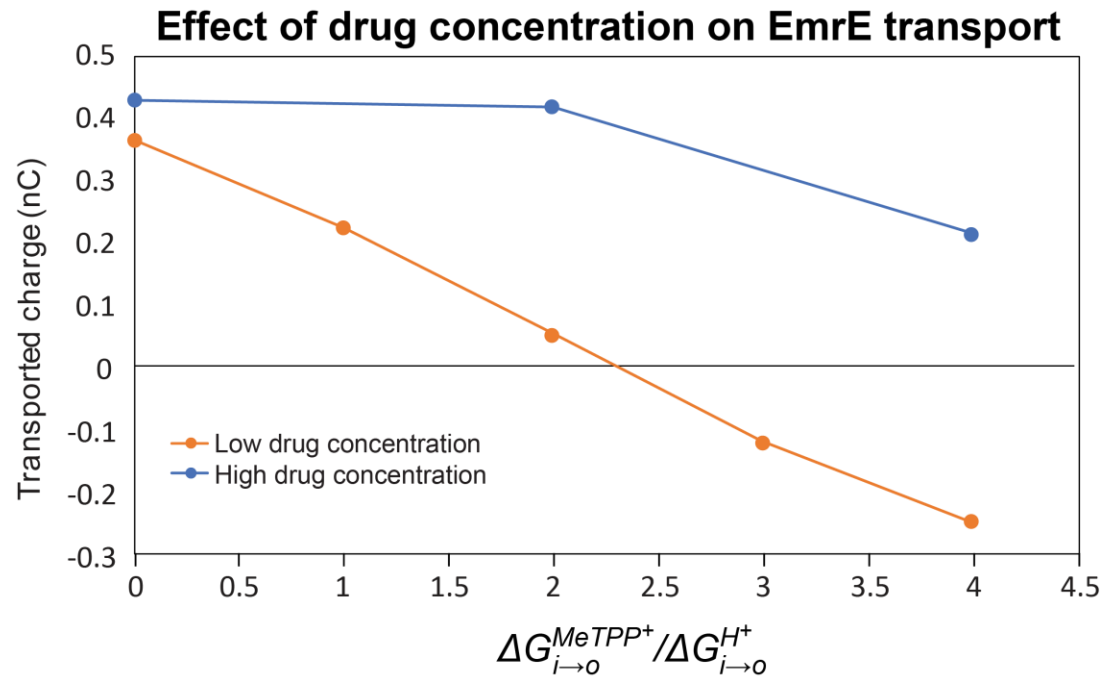
Sensor internal solution can be changed

$$R^2 = 0.999$$

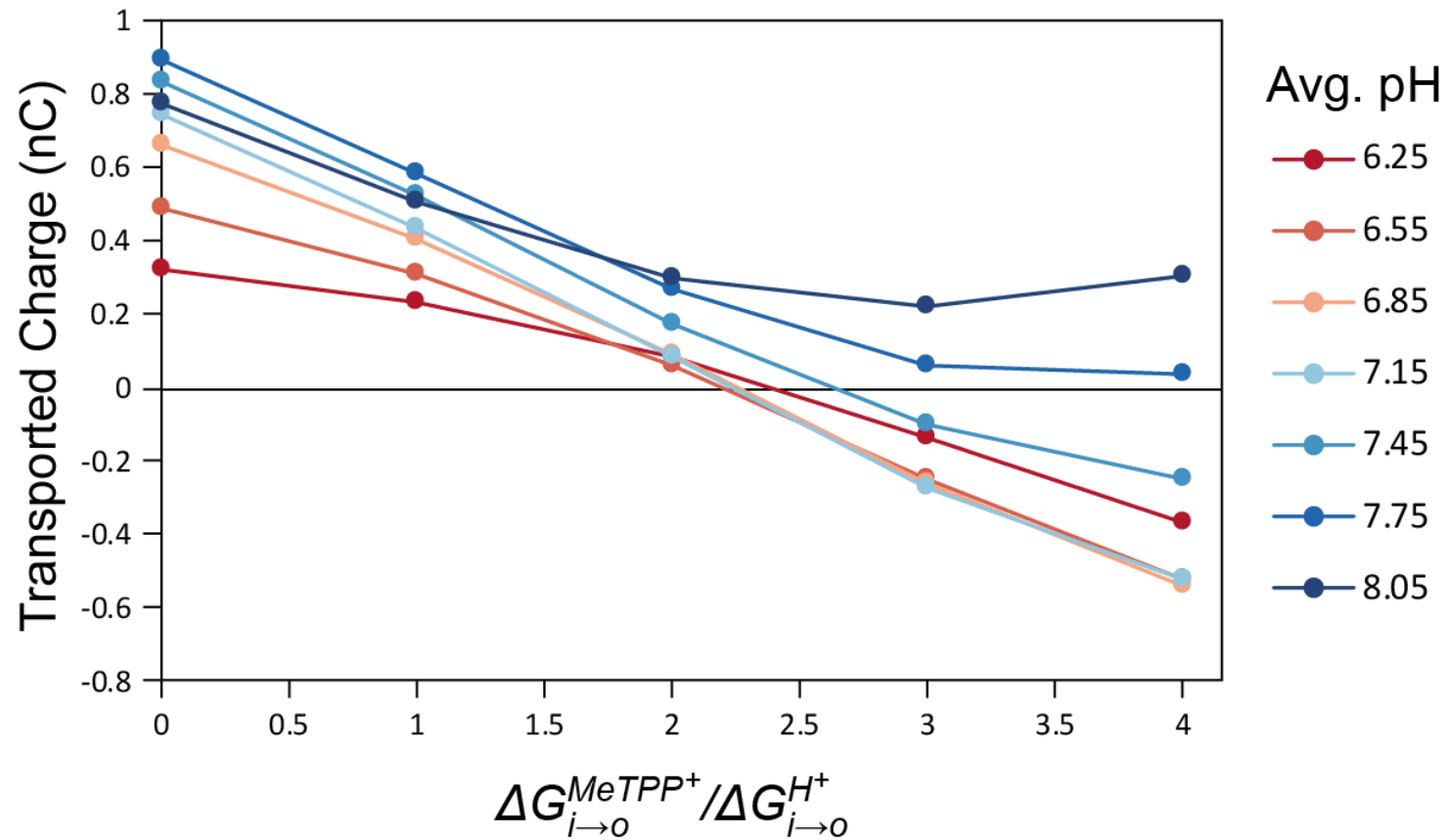
$$\text{x-intercept} = 2.02 \pm 0.06$$



EmrE has variable stoichiometry



EmrE has variable stoichiometry



35 experimental conditions tested <3 hours, using <50 picomol protein

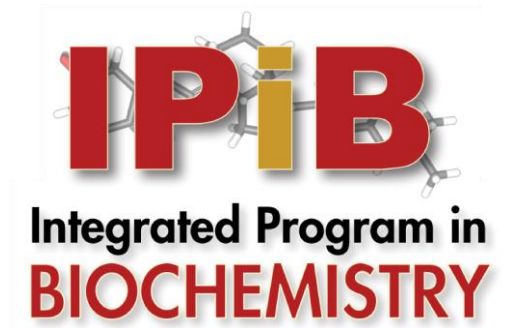
Advantages of SURFE²R Reversal Potential

- No need for fluorescent probes or radioactive isotopes
- Much higher throughput – hundreds of measurements in a day
- Requires very little sample
 - 10^{-10} mol protein per sensor
 - Up to 100 measurements per sensor
 - Can measure multiple conditions on each sensor

Acknowledgements

Henzler-Wildman Lab

Grant Hisao
Vilius Karauskas
Adam Lewis
Peyton Sprecker
Andrea Killian



Nanion Technologies

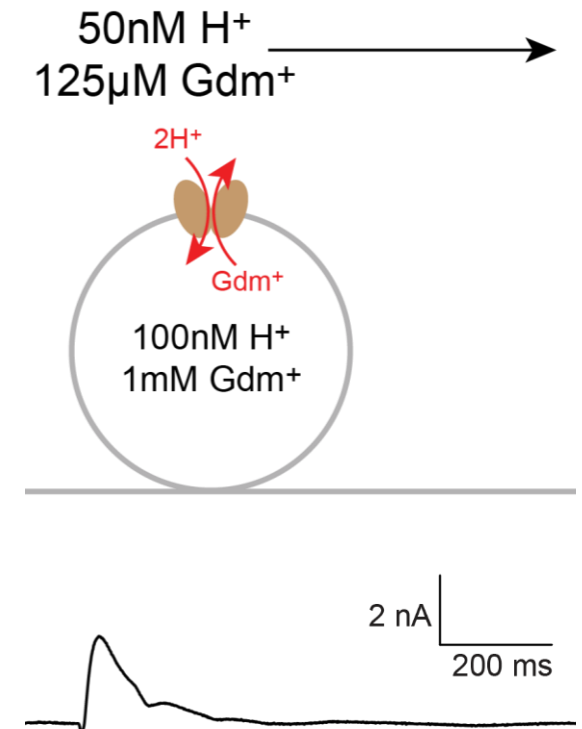
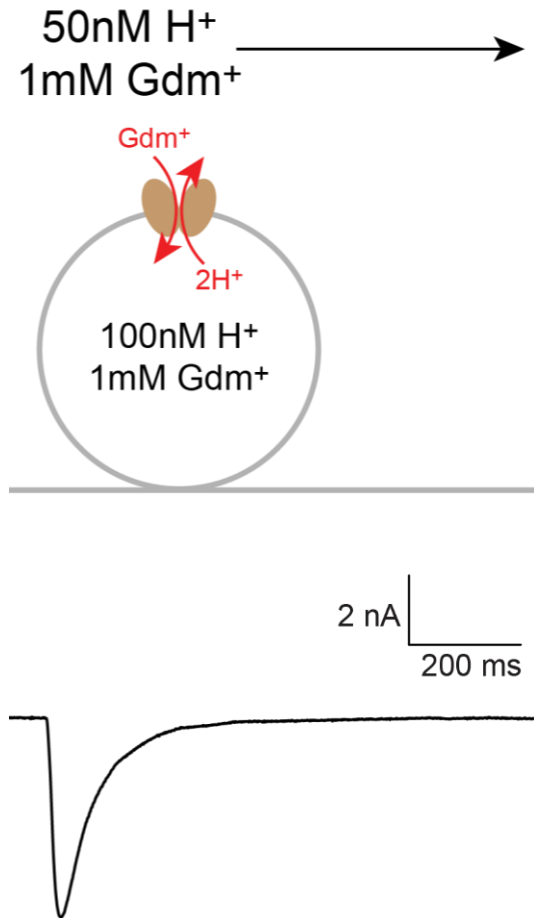
Dr. Maria Barthmes
Dr. Andre Bazzone



Funding

NIH 1R01GM095839 and T32 GM007215
Nanion SURFE²R N1 Grant 2018

SURFE²R Reversal Potential



Derivation of reversal potential equation

For coupled antiport:

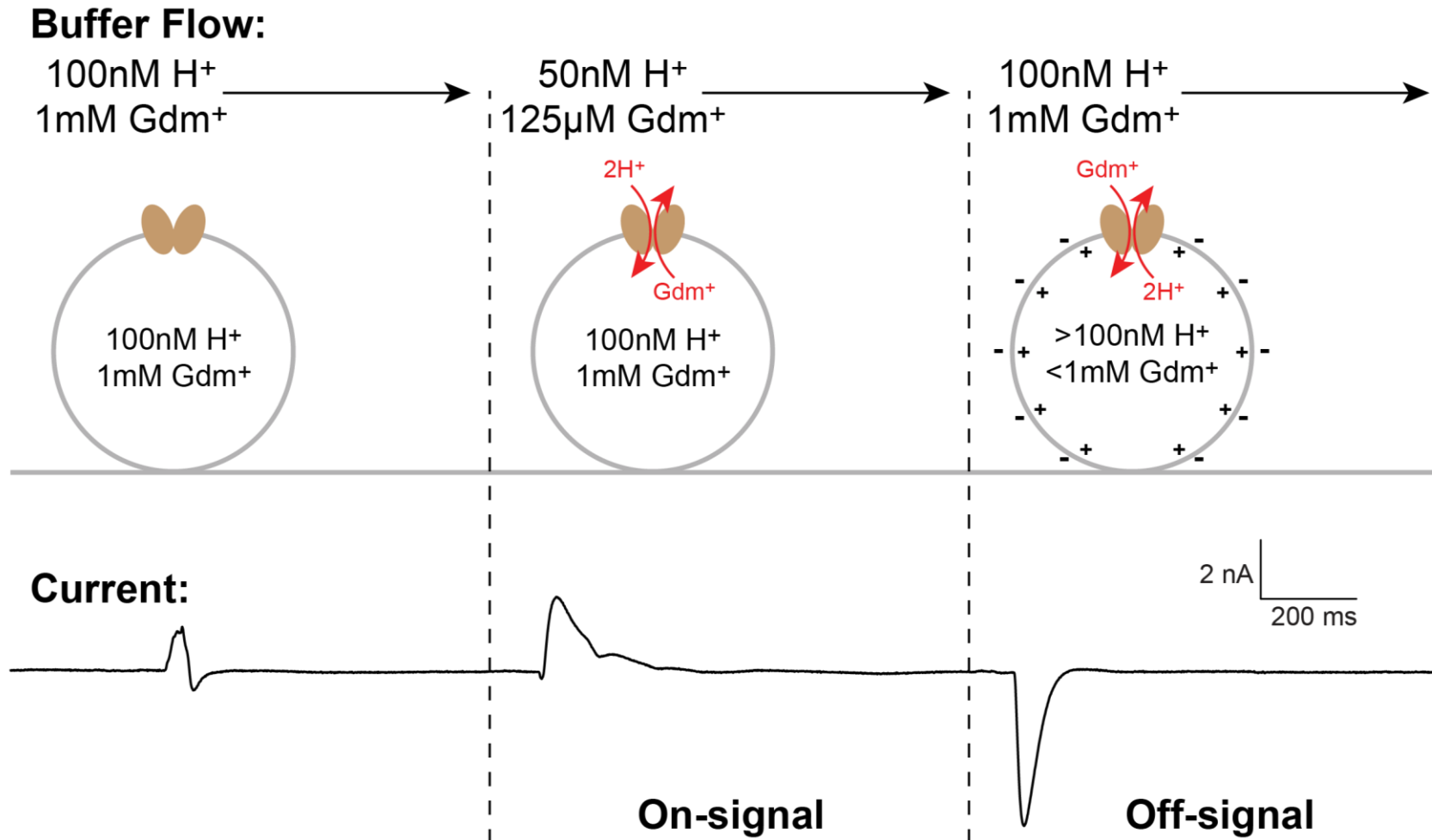


Eq. 2 $\Delta G_{\text{antiport}} = n\Delta G_{i \rightarrow o}^{\text{Gdm}^+} - m\Delta G_{i \rightarrow o}^{\text{H}^+}$

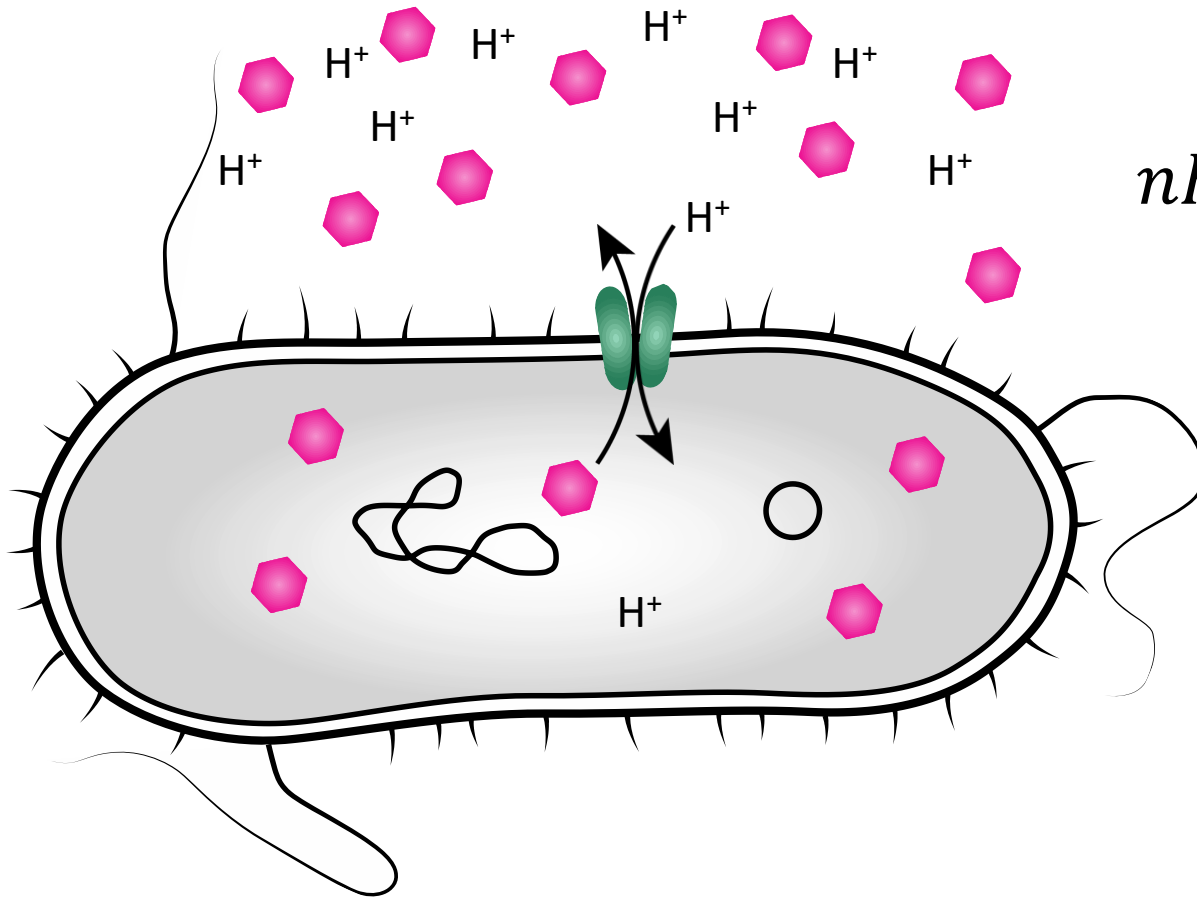
When $\Delta G_{\text{antiport}} = 0$,

Eq. 3 $\frac{m}{n} = \frac{\Delta G_{i \rightarrow o}^{\text{Gdm}^+}}{\Delta G_{i \rightarrow o}^{\text{H}^+}}$

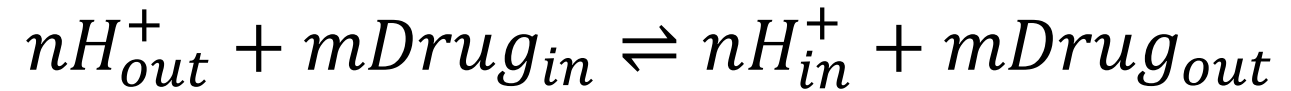
Assay Scheme – Guanidinium gradient drives proton transport



Thermodynamics of proton coupled drug efflux



Antiport



$$\left(\frac{H_{out}^{+}}{H_{in}^{+}} \right)^n = \left(\frac{Drug_{out}}{Drug_{in}} \right)^m$$

***E. coli* EmrE is a model drug/proton antiporter**

SSM Reversal Assay is quantitative

