

Ion Channel Drug Discovery – Recent Advances in Novel Non-Opioid Pain Research

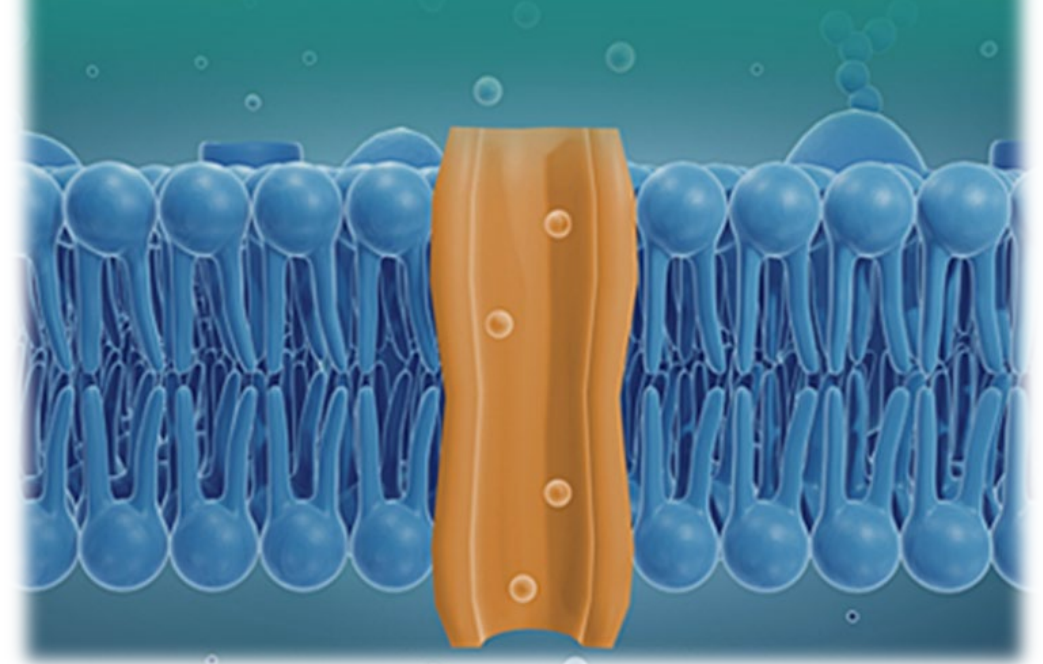
John Gilchrist, Ph.D., Principal Scientist
Latigo Biotherapeutics Inc.



James Costantin, Ph.D., Scientific Market Development Manager – Ion Channels
Eurofins DiscoverX®

1. Ion Channels in Drug Discovery and Development
 - The Pain Pathway and Current Therapeutics
2. Case Study – “LTGO-33 Is a Novel Nav1.8 Inhibitor”
 - Presented by John Gilchrist – Latigo Biotherapeutics Inc.
3. Eurofins DiscoverX[®] Ion Channel Products
 - Capabilities, Stable Cell Lines, and Ready-to-assay Cells
4. Summary and Conclusions

Ion Channel Drug Discovery – Recent Advances in Novel Non-Opioid Pain Research



Eurofins DiscoverX[®] a Global Leader in Cell-based Assays for Screening, and Profiling



DiscoverX

Accelerating Breakthroughs from Discovery to Delivery

25+ Years of Experience Building Cell-based Assays



Fremont, California | Poitiers, France | Shanghai, China



CALIXAR

Pure & Native Full Length
Membrane Proteins

Functional Kinases

GPCRs

Industry's Largest Cell-based Assay Portfolio

10+ Druggable Target Classes

60+ Ion Channel Assays

Stable Cell Lines

Ready-to-Assay

Membrane Preps

Target Discovery

Hit Screening

Lead Optimization

Safety Studies

Internally Validated

>30 Billion

Data Points

Ion channels represent a key target class for therapeutics to treat human disease

- Over 400 genes encode for ion channels and their accessory subunits in humans
- They are ubiquitous proteins found in all cell types
- Ion channels are involved in normal physiological processes such as the transmission of pain signals in the central and peripheral nervous system
- Therapeutic intervention in normal or aberrant ion channel function can lead to effective treatment of human diseases, including the treatment of pain disorders

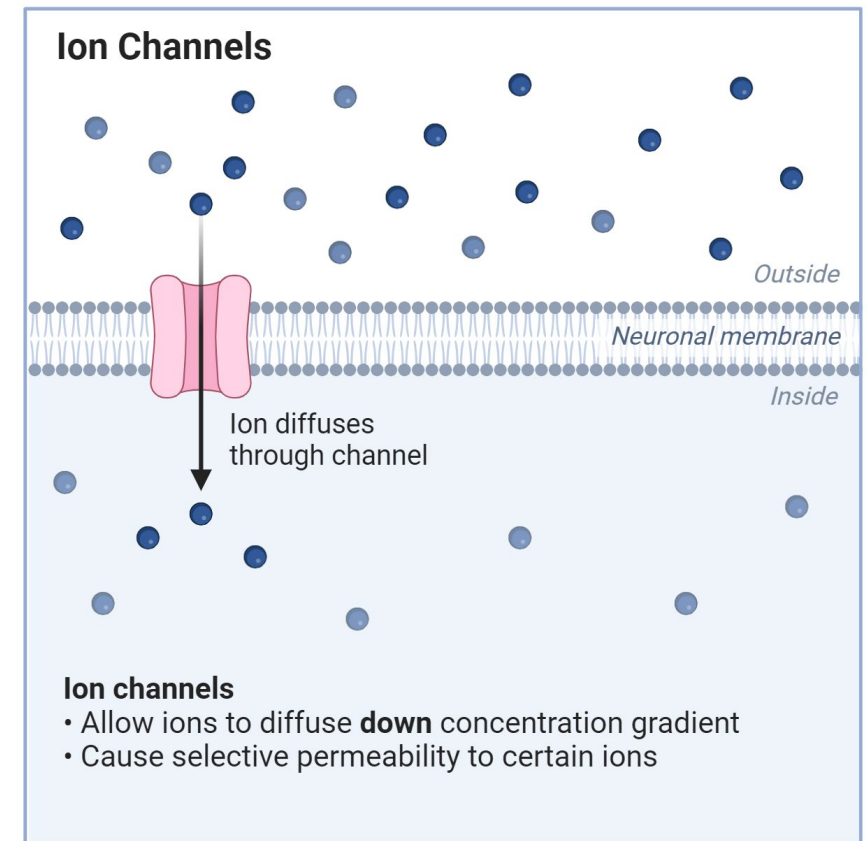


Image source: BioRender.com

Channelopathies

- Mutations in ion channel genes or their related proteins cause diseases (channelopathies) that range from mild to severe, affecting the nervous, cardiovascular and respiratory systems, as well as causing endocrine, kidney and immune-related diseases

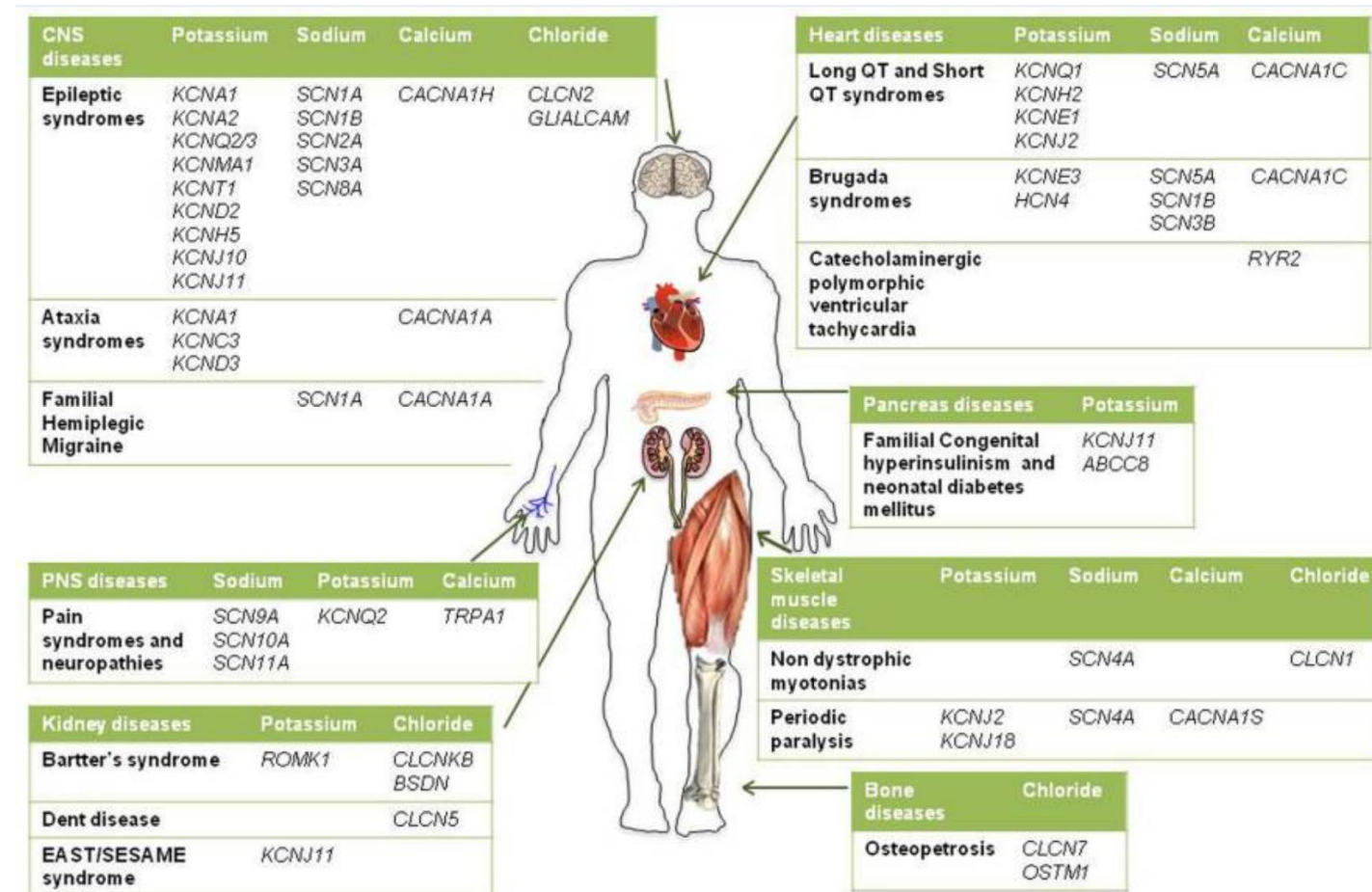


Image source: Front. Pharmacol., 09 May 2016, Vol.7.
<https://doi.org/10.3389/fphar.2016.00121>

The pain pathway, opioids, and the need for non-addictive analgesics

The Pain Pathway

- Pain stimuli are detected by nociceptors in the peripheral nervous system
- Signals are then transmitted to the Central Nervous System (CNS) via Dorsal Root Ganglion (DRG) neurons
- Sodium channels in DRG neurons initiate trains of action potentials in response to the incoming peripheral nerve signals
- This transmits the pain signal to the brain where it is perceived

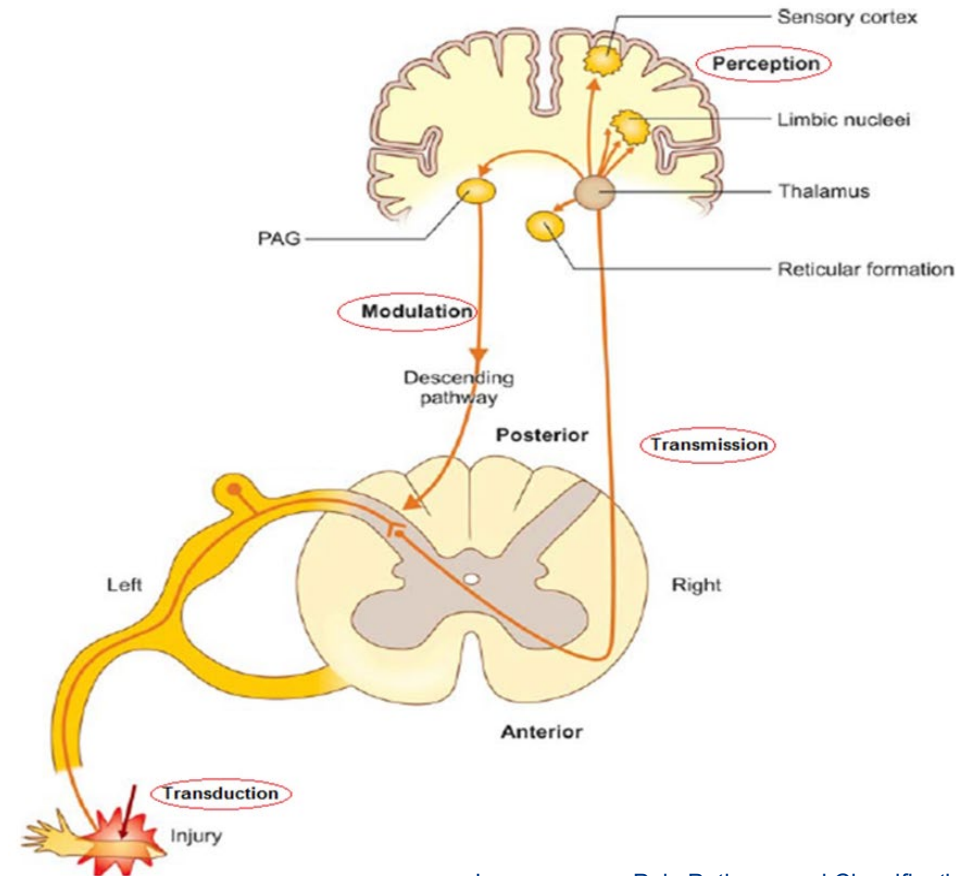
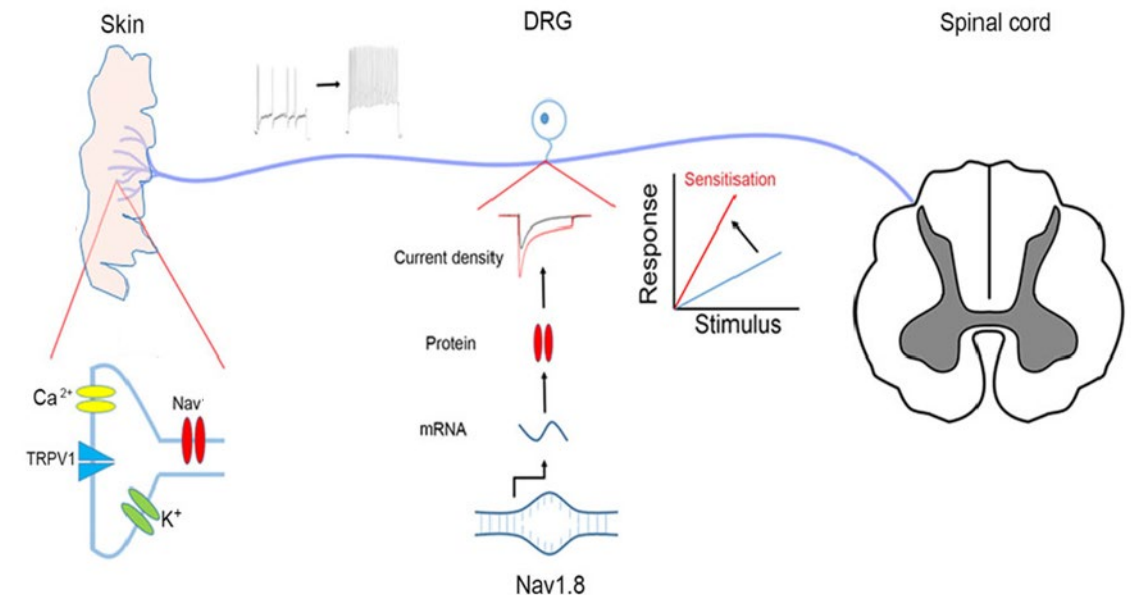


Image source: Pain Pathway and Classification of Pain (Daradia, The Pain Clinic <https://daradia.com/pain-pathway/>)

The pain pathway, opioids, and the need for non-addictive analgesics

- Inhibition of Nav1.8 ion channels in DRG neurons should be an effective analgesic by reducing transmission of the pain signal from DRG neurons to the brain
- Highly addictive opioid drugs act by depressing neuronal excitability in the CNS by reducing the activity of Ca^{2+} channels and increasing the activity of K^{+} channels
- Development of an effective inhibitor of Nav1.8 ion channels could lead to a non-addictive analgesic



Adapted from: *Acta Biochim Biophys Sin*, 2016, 48(2), 132–144. doi: 10.1093/abbs/gmv123



LATIGO BIO

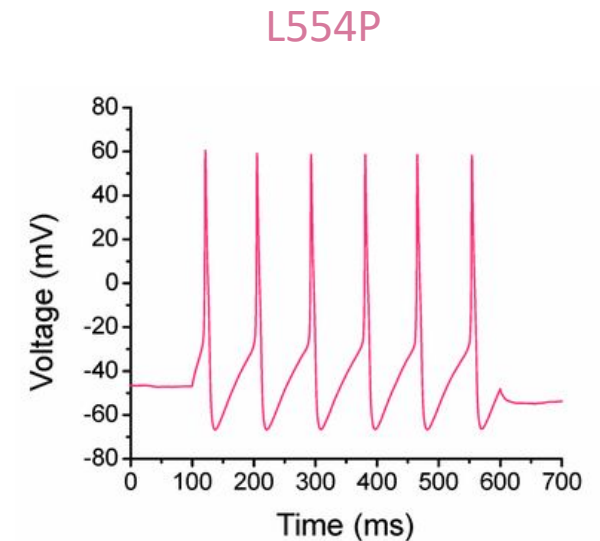
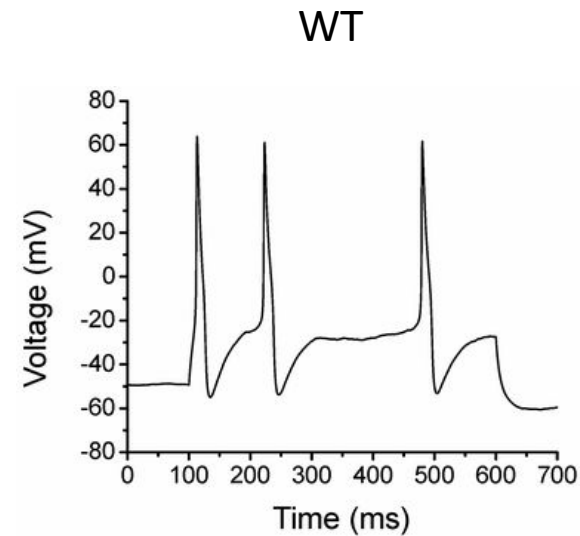
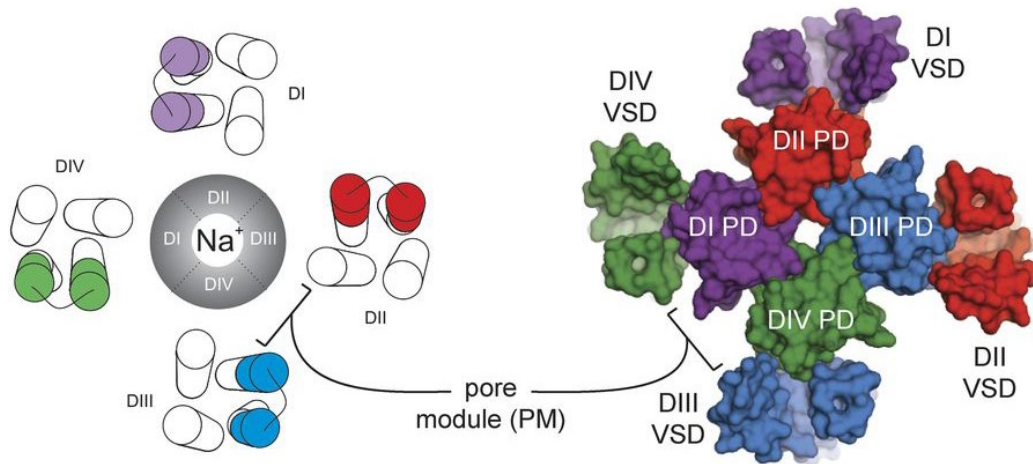
LTGO-33 Is a Novel $\text{Na}_v1.8$ Inhibitor

John Gilchrist, Ph.D.

Latigo Biotherapeutics Inc.

15 May 2024

Na_v1.8 Is a Validated Target for Analgesia



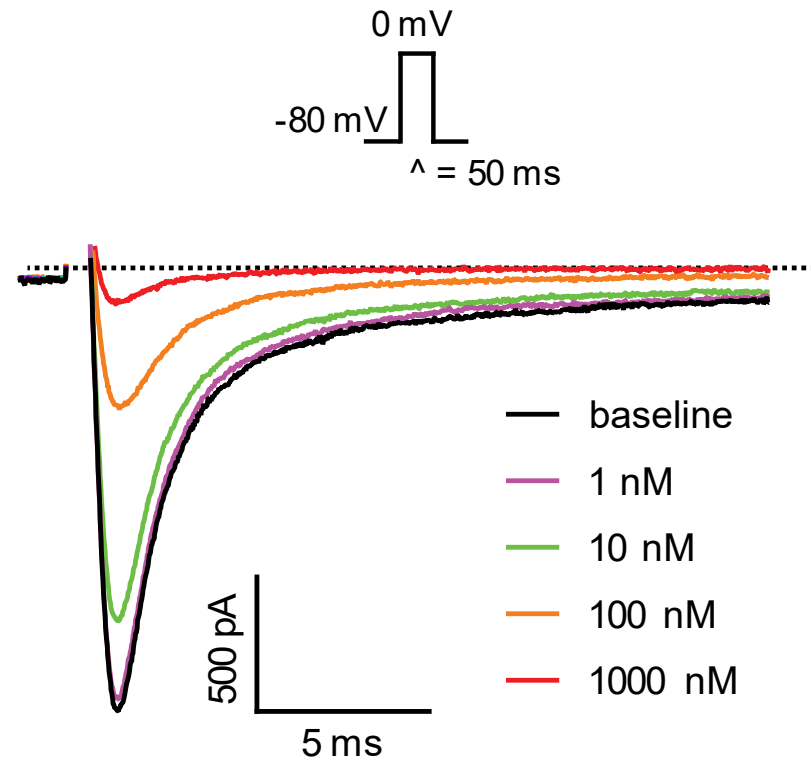
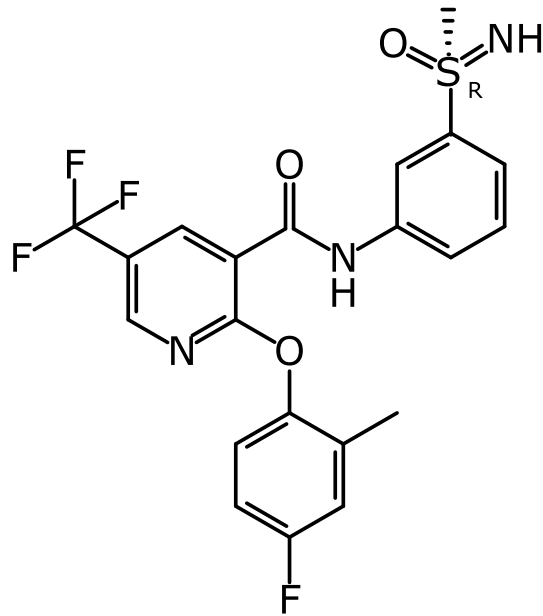
Na_v1.8 is a voltage-gated sodium channel found in nociceptive DRG sensory neurons

Genetically- and pharmacologically-validated as a target for pain

Ahern, Christopher A., et al. "The hitchhiker's guide to the voltage-gated sodium channel galaxy." *Journal of General Physiology* 147.1 (2016): 1-24.

Faber, Catharina G., et al. "Gain-of-function Nav1.8 mutations in painful neuropathy." *Proceedings of the National Academy of Sciences* 109.47 (2012): 19444-19449.

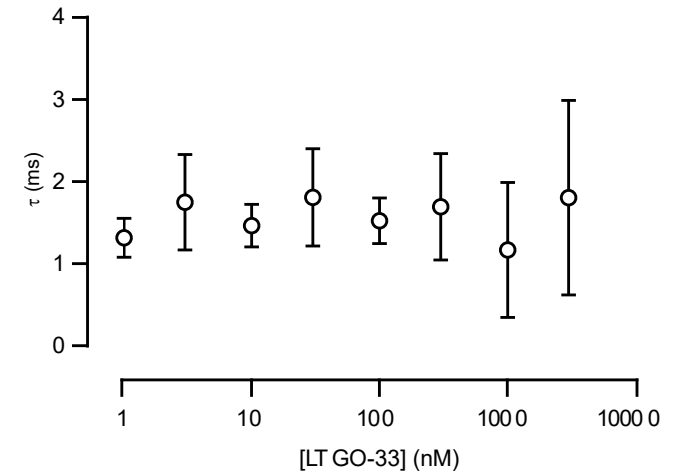
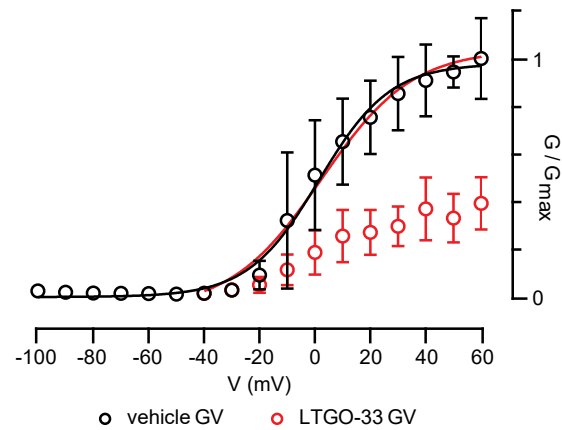
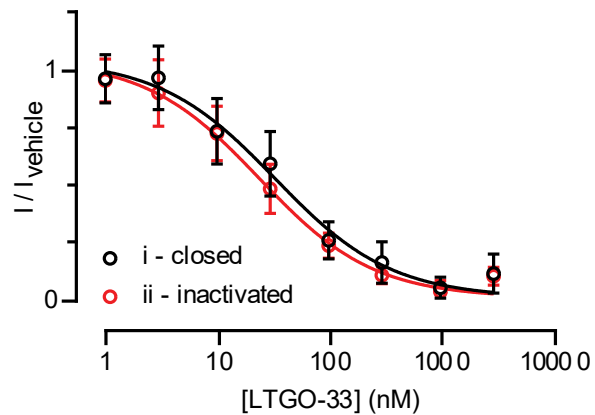
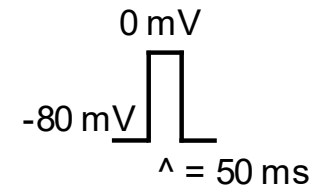
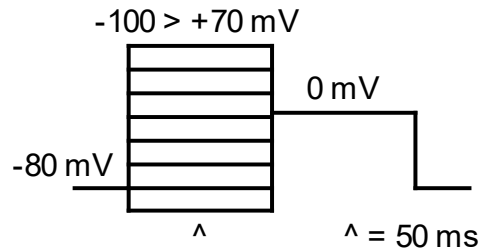
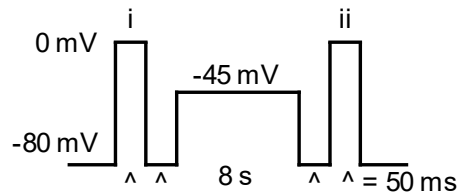
LTGO-33 Is a Potent $\text{Na}_v1.8$ Inhibitor



LATIGO BIO

LTGO-33 inhibits $\text{Na}_v1.8$ with nM potency

LTGO-33 Inhibition Is State-Independent

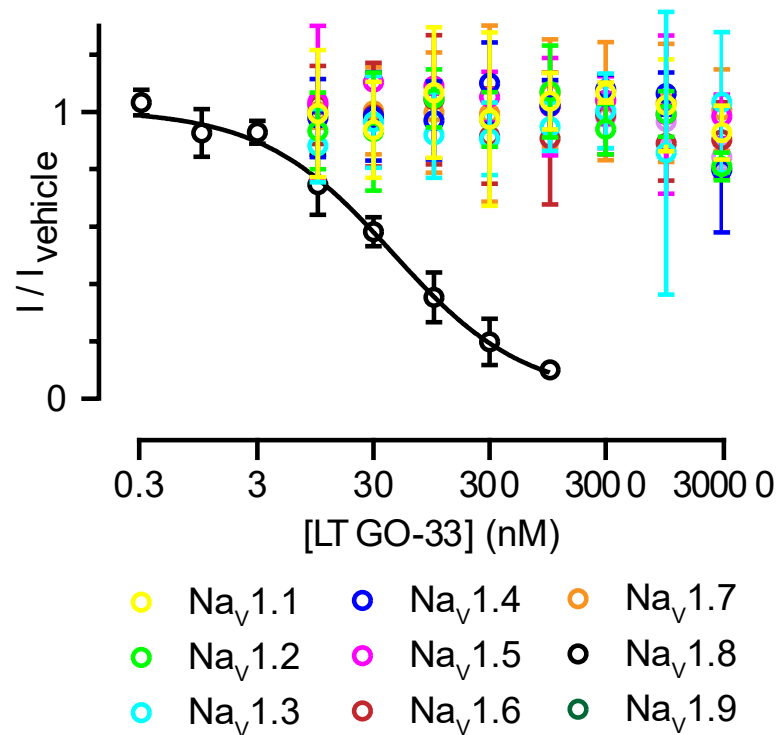


No preference for inhibition of resting vs inactivated state

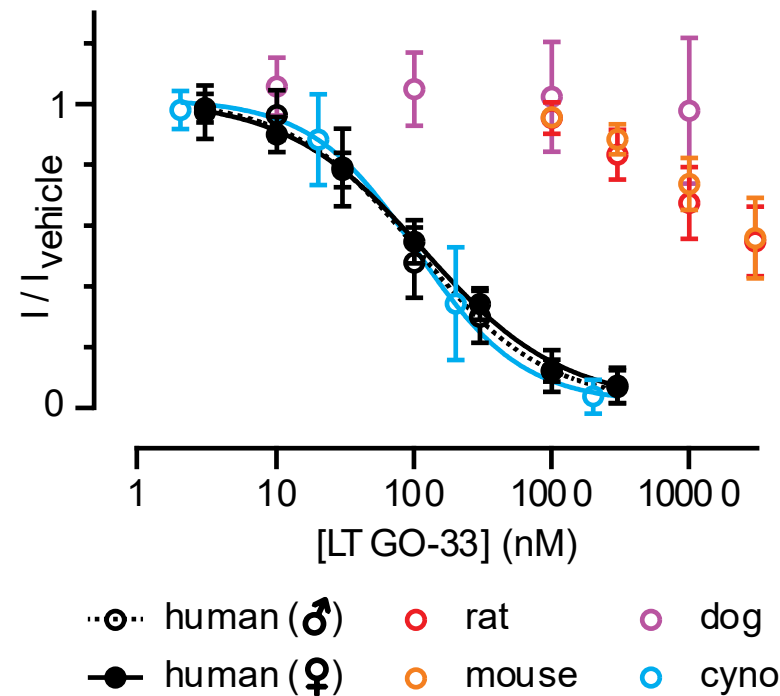
No shift in voltage-dependence of activation

No change in rate of fast inactivation

LTGO-33 Is Isoform- and Species-selective

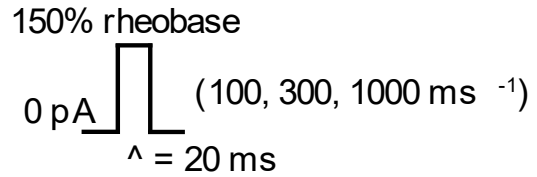


>600-fold selective for Na_v1.8 vs Na_v1.1-1.7 and Na_v1.9

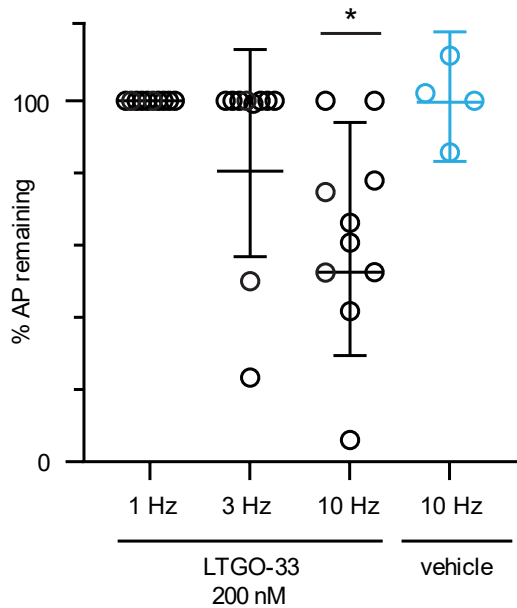


>250-fold selective for primate TTX-R vs rodent and dog TTX-R

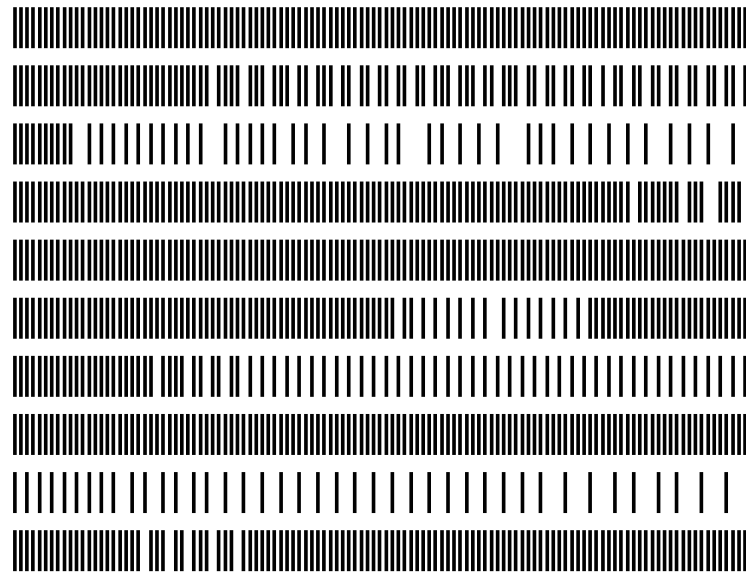
LTGO-33 Inhibits Action Potentials in hDRG Neurons



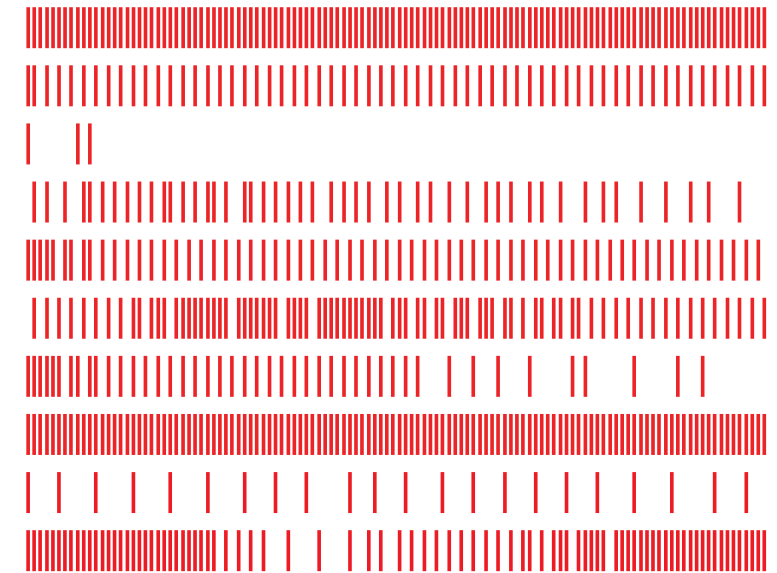
TTX-free to simulate physiological conditions



baseline (10 Hz)

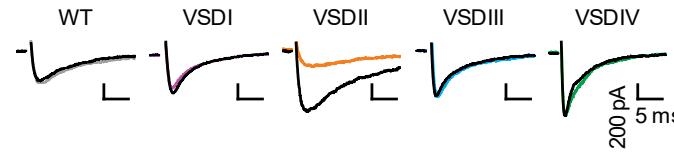
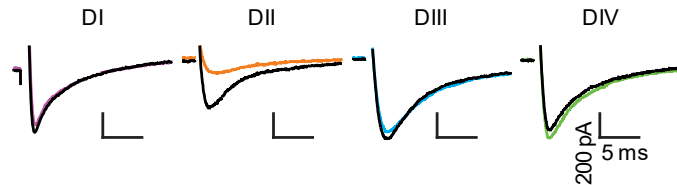
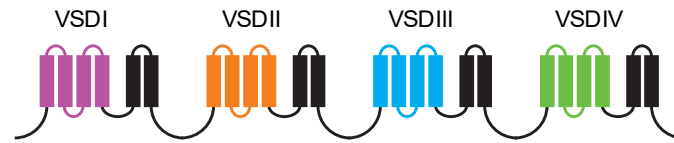
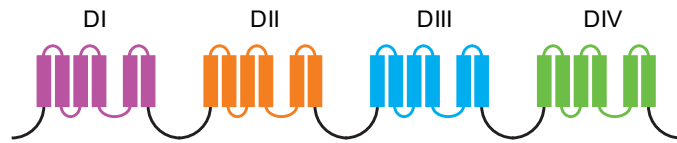


200 nM LTGO-33 (10 Hz)



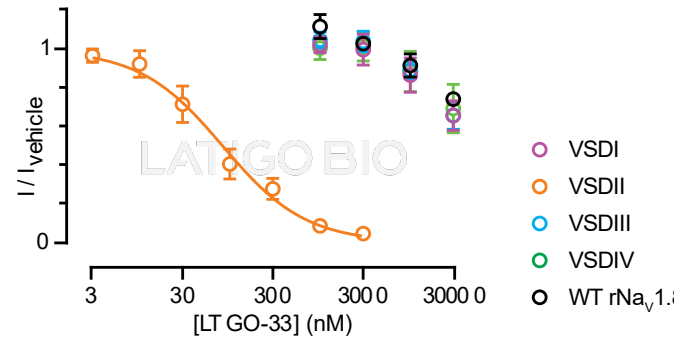
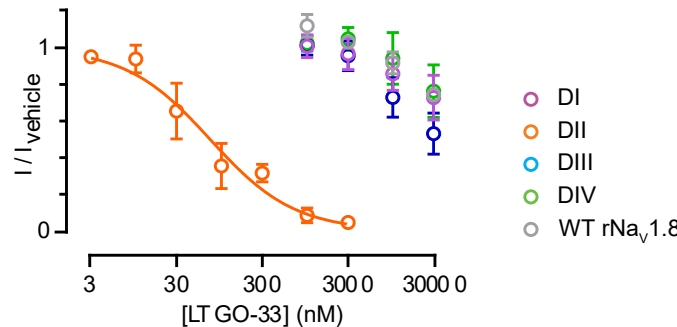
LTGO-33 is effective at decreasing AP firing at a higher frequency (10 Hz) at 2x hDRG IC₅₀

VSDII Is Required for LTGO-33 Activity



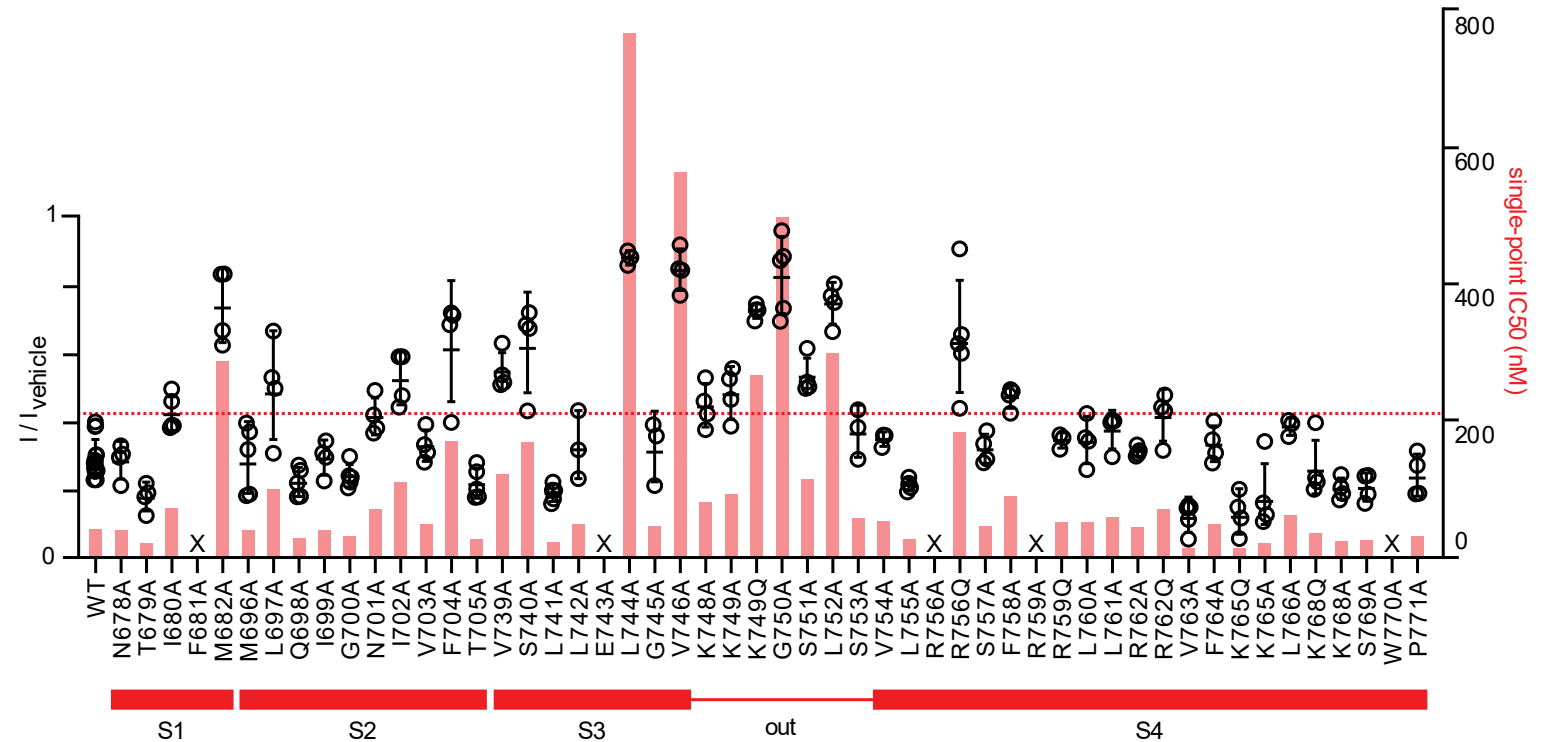
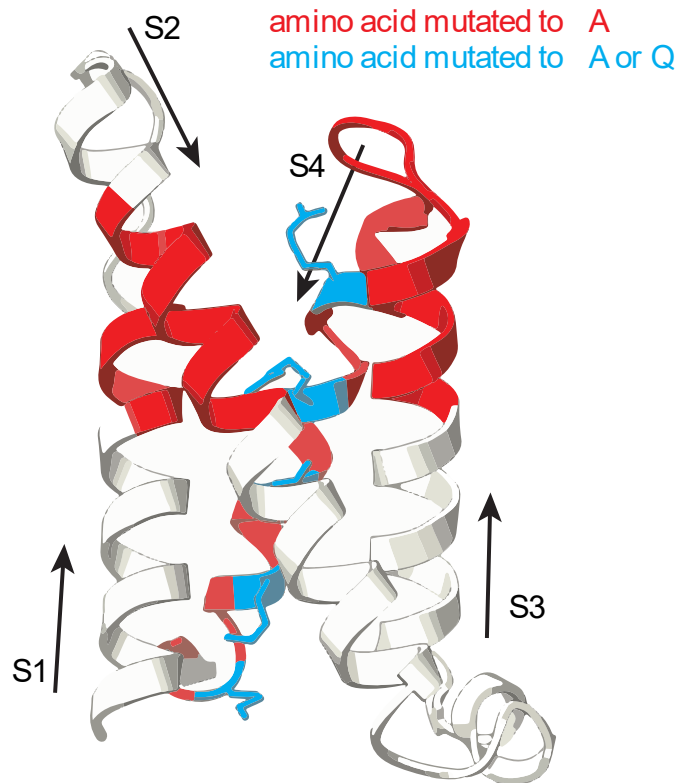
rNa_v1.8 (with reduced LTGO-33 sensitivity) used as host for gain-of-function domain chimeras

DII (IC₅₀ 74 nM) and VSDII (IC₅₀ 84 nM) chimeras inhibited by LTGO-33, all others IC₅₀ >30 μM



Comparable potency on DII and VSDII chimeras suggests VSD sufficient, pore unnecessary

Residues in S1, S3, and S4 Critical for LTGO-33 Activity

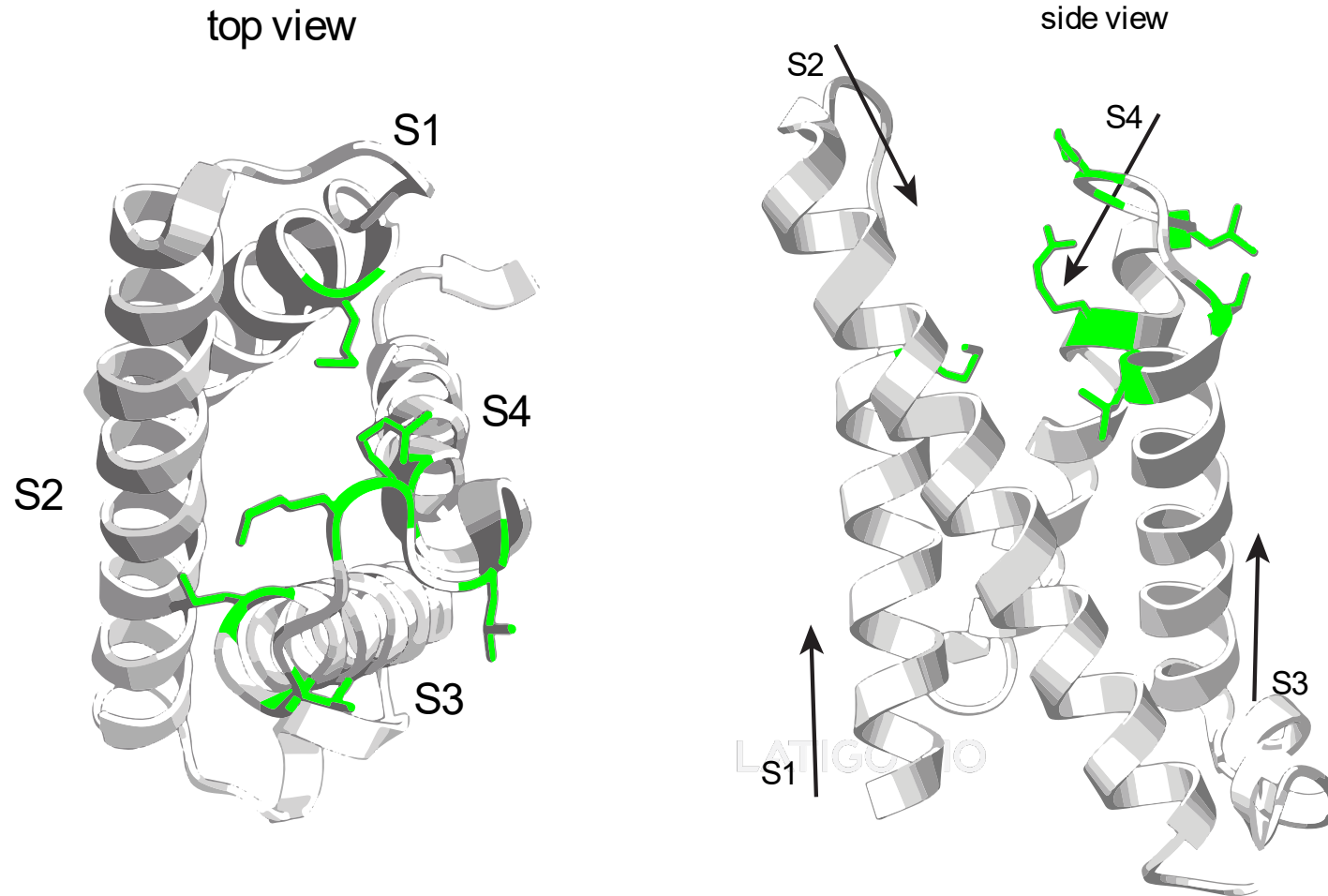


LATIGO BIO

Alanine scan conducted through regions of dissimilarity between human and rat

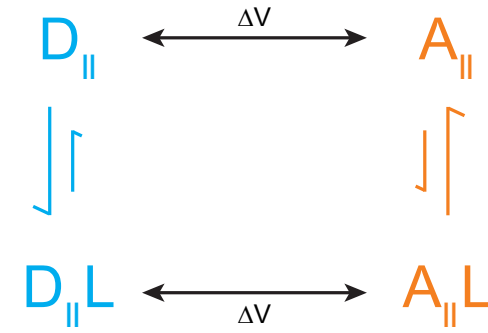
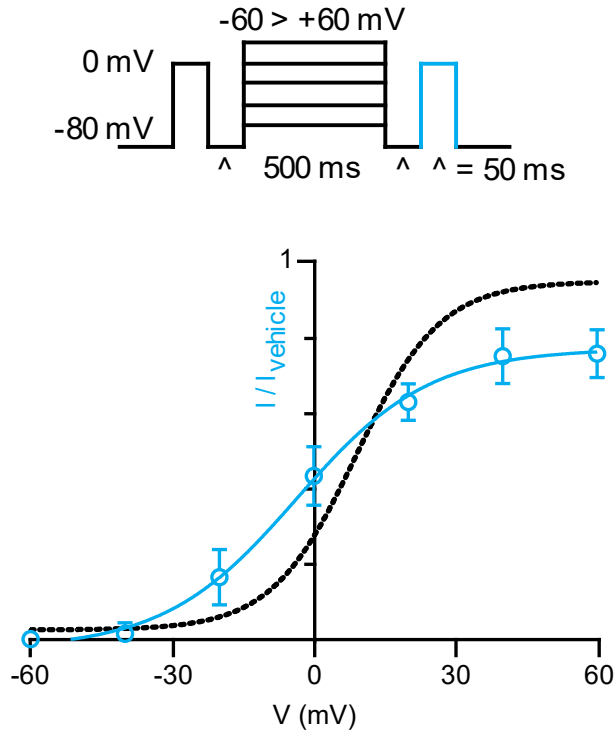
Amino acids showing >5x decrease in potency found in S1 and S3b-S4 loop

LTGO-33 Interaction Site in Extracellular VSDII



Important residues found in outer portion of VSDII – similar to ProTx-II binding site

Disinhibition and a Gating Model

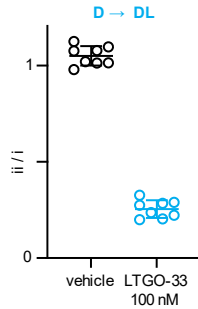
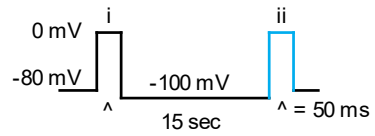


A basic model proposes LTGO-33 stabilizes the VSDII deactivated (down) state to prevent channel activation

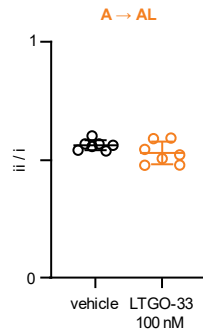
Depolarization can relieve Na_v1.8 inhibition by LTGO-33, akin to ProTx-II

LATIGO BIO

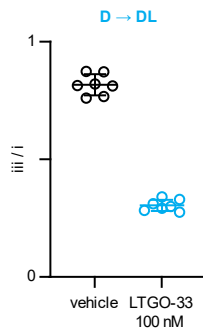
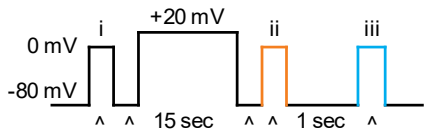
LTGO-33 Inhibits Closed but not Activated Channels



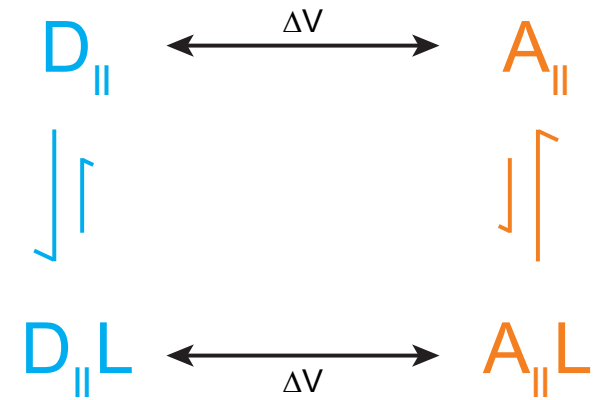
LTGO-33 can inhibit channels from the closed state (held at negative Vm)



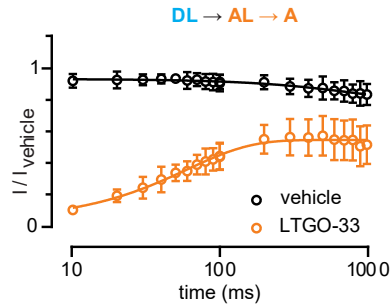
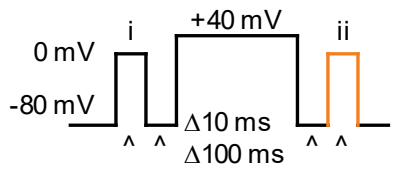
Similar inhibition in vehicle- and LTGO-33-treated cells due to slow inactivation, minimal further inhibition with LTGO-33



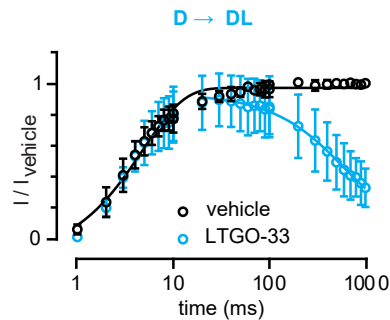
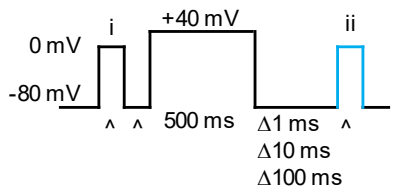
LTGO-33 inhibits channels after brief hold at negative Vm (closed state)



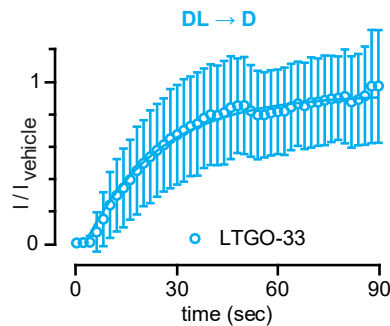
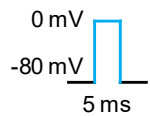
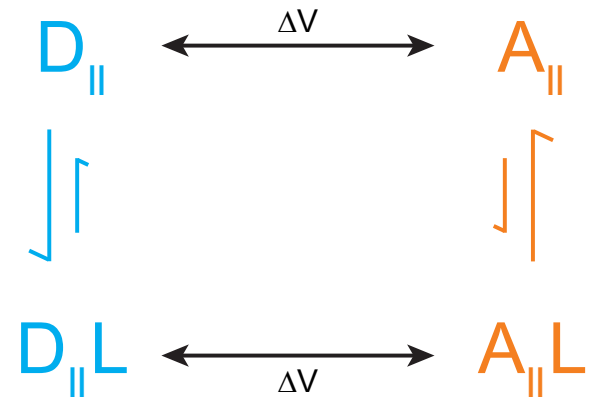
LTGO-33 Inhibition Can Be Relieved by Depolarization



LTGO-33 disinhibition is complete within a 1 second depolarization

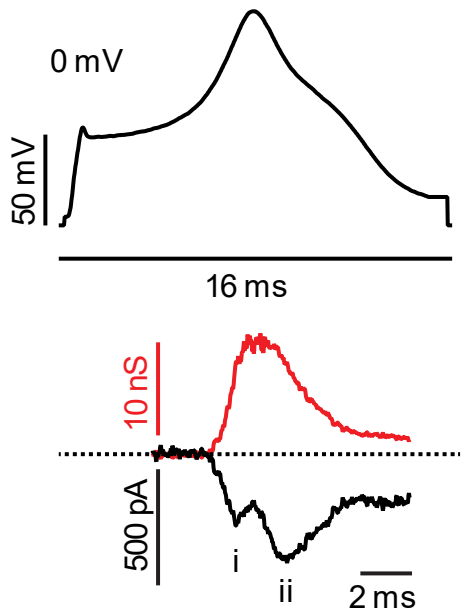


LTGO-33 reinhibition occurs rapidly in the deactivated state

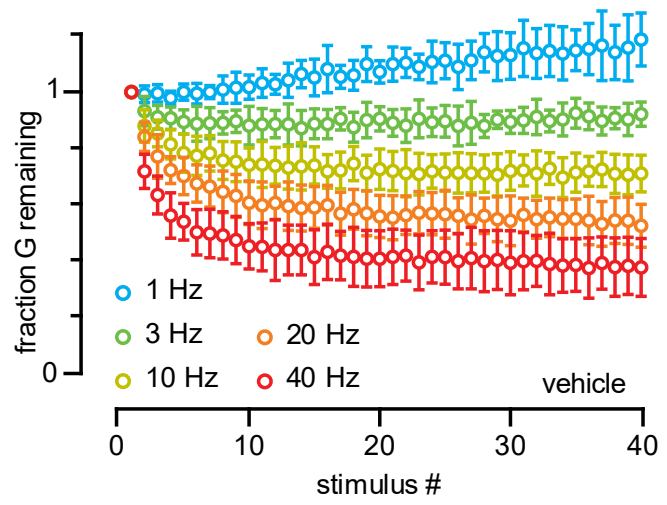


LTGO-33 is relatively slower to dissociate from deactivated VSDII

LTGO-33 Exhibits Minimal Reverse-use Dependence

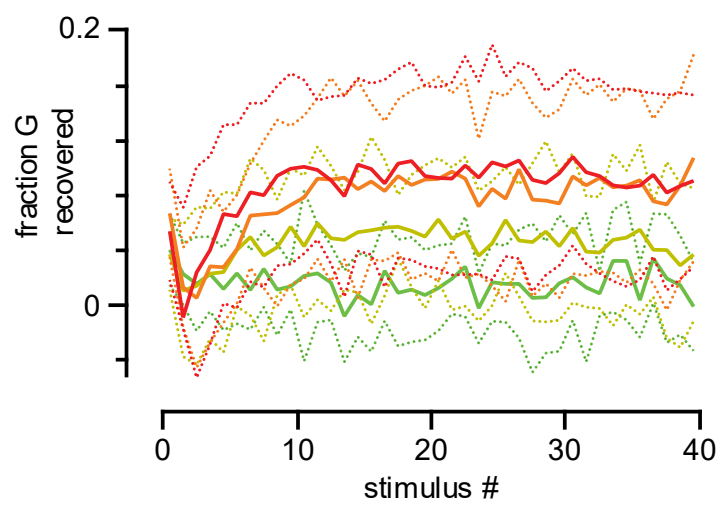
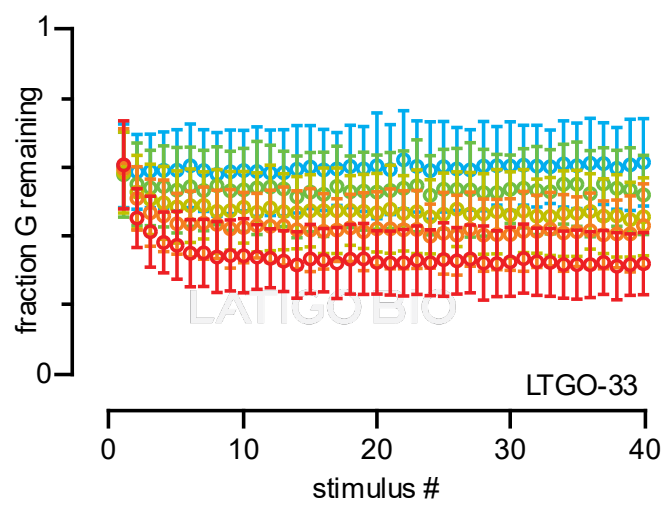


AP waveform applied to ND7/23 cells transfected with hNa_v1.8, conductance plotted

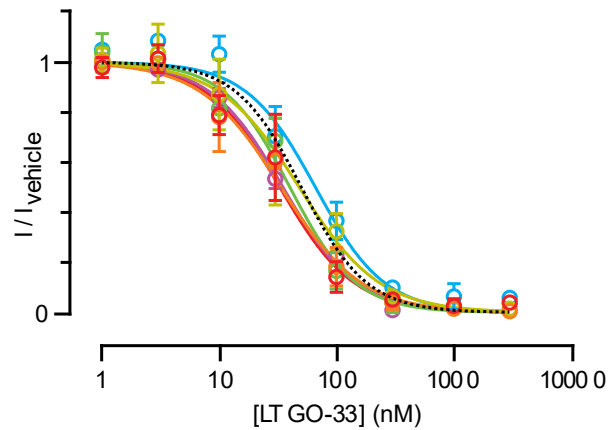
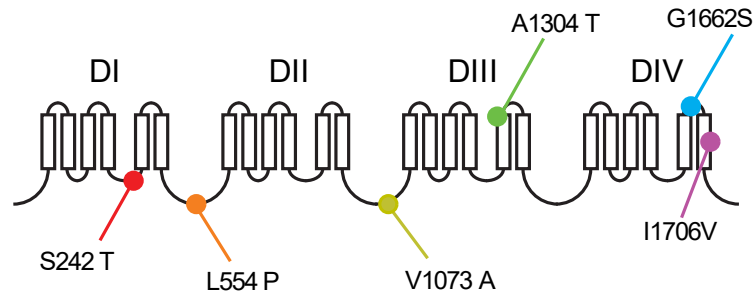


Use-dependent inhibition (UDI) observed at higher frequencies

Curve-fitting approach subtracted UDI from LTGO-33 treated cells – minimal reverse-use dependence observed up to 40 Hz



LTGO-33 Inhibits Nav1.8 Harboring Variants Associated with Pain Disorders



- S242T
- L554P
- V1073A
- A1304T
- G1662S
- I1706V

LATIGOBIO

WT hNav_v1.8

Na_v1.8 variants associated with pain disorders located outside of VSDII do not affect LTGO-33 pharmacology

Conclusions

- LTGO-33 is a potent, selective Na_v1.8 inhibitor that blocks AP firing in an *in vitro* human DRG neuronal model
- Unlike prior published Na_v1.8 inhibitors, LTGO-33 shows no preference for inhibiting from inactivated state
- LTGO-33 engages the channel at a novel site in VSDII similar to peptide spider toxins
- LTGO-33 is effective in inhibiting Na_v1.8 variants associated with pain disorders

Acknowledgements

Latigo Team

Victoria Jiang
Nien-Du Yang
Bryan Moyer

Shanti Amagasu
Tina Holt

External Activities

AnaBios
Neuroservice USA
Charles River Labs
B'SYS
GenScript

Find the paper at *Molecular Pharmacology!*

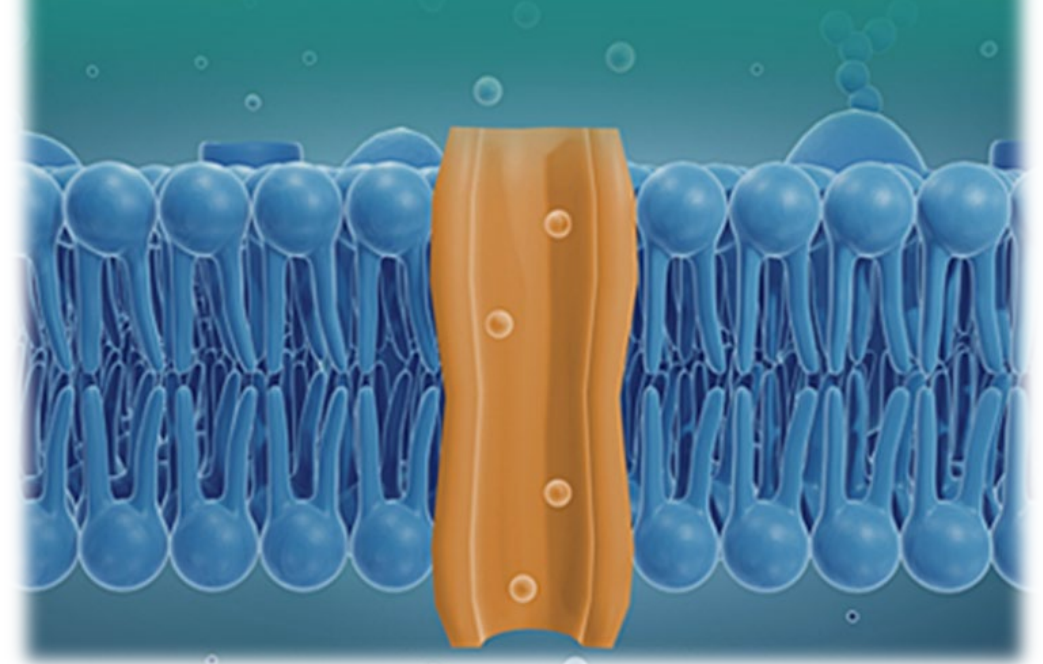
The hNa_v1.8/β1 (CYL3025) and rNa_v1.8 (CYL3050) stable cell lines are products of Eurofins DiscoverX.

LATIGO BIO



1. Ion Channels in Drug Discovery and Development
 - The Pain Pathway and Current Therapeutics
2. Case Study – “LTGO-33 Is a Novel Nav1.8 Inhibitor”
 - Presented by John Gilchrist – Latigo Biotherapeutics Inc.
3. Eurofins DiscoverX[®] Ion Channel Products
 - Capabilities, Stable Cell Lines, and Ready-to-assay Cells
4. Summary and Conclusions

Ion Channel Drug Discovery – Recent Advances in Novel Non-Opioid Pain Research



Accelerating Breakthroughs from Discovery to Delivery

Ion Channel Stable Cell Lines Used for Drug Discovery for 15+ Years



Fremont, California | Lyon, France



CALIXAR

Functional Kinases,
GPCRs and Ion
Channels

Reconstituted Full Length
Membrane and Soluble
Proteins

Nanodiscs, Proteoliposomes,
and Amphipols

65+ Ion Channel
Assays

Stable Cell Lines

Ready-to-Assay Frozen
Cells

Membrane Preps

Membrane Trafficking
Assays

Target Discovery

Hit Screening

Lead Optimization

Safety Studies

Validation Data by
Electrophysiology
Available for Cell
Lines

PrecisION™ Ion Channel Cell Lines

- High-quality cell lines for target discovery, hit screening, lead optimization, and safety studies
- Suitable for manual and automated electrophysiology and fluorescence imaging studies
- Available as continuous culture and now in a ready-to-assay (RTA) format

Eurofins DiscoverX is now the home and exclusive provider of PrecisION Ion Channel cell lines

- Eurofins DiscoverX, part of Eurofins Discovery, the products company and the leading provider of cell lines, cell-based assays, and enzymes for drug discovery and development
- PrecisION ion channel stable cell lines portfolio was acquired through the acquisition of Millipore (EMD Millipore/Merck Millipore) after a series of smaller acquisitions

Cell line production, marketing, and technical support

- Quality control (QC) validation of existing Eurofins DiscoverX PrecisiON™ cell lines at the Fremont, CA lab facility
- Development and QC validation of frozen RTA PrecisiON channel cell lines
- Validation of custom ion channel targets through custom development capabilities program
- Expansion of our Custom Development Capabilities program into rare disease targets
 - Channelopathy variants and wild type controls



Nanion SyncroPatch 384i

PrecisION[™] channel cell lines

hERG (K⁺)

hERG-CHO
hERG-HEK

K⁺ Channels

K_v1.1-CHO
K_v1.2-CHO
K_v1.3-CHO
K_v1.4-CHO
K_v1.5-CHO
K_v1.6-CHO
K_v1.7-CHO
K_v1.8-CHO
K_v12.2-HEK
K_v2.1-CHO
K_v2.1/K_v9.2-CHO
K_v3.1-CHO
K_v3.2-CHO
K_v3.3-CHO
K_v4.2/KChIP2-CHO

K⁺ Channels

K_v4.3/KChIP1-CHO
K_v4.3/KChIP2-HEK
K_v7.2/K_v7.3-CHO
K_v7.3/K_v7.5-CHO
K_v7.4-HEK
K_v7.4/K_v7.5-HEK
K_{ir}2.1-HEK
K_{ir}6.2/SUR2A-HEK
KNCQ1/hminK-CHO

HCN Channels

HCN1-HEK
HCN2-HEK
HCN3-HEK
HCN4-CHO

Ca²⁺ Channels

Ca_v1.2 $\alpha_{1C}/\beta_{2a}/\alpha_2\delta_1$
Ca_v2.2-HEK
Ca_v3.2-HEK

Other Channels

ASIC3-HEK
CFTR-HEK
TRPA1-HEK
TRPV1-HEK
TRPV3-HEK
TRPV4-HEK

Ligand-gated Channels

nAChR $\alpha_1/\beta_1/\delta/\epsilon$ -HEK
nAChR α_3/β_4 -HEK
nAChR α_4/β_2 -HEK
nAChR $\alpha_7/ric3$ -HEK
GABA_A $\alpha_1/\beta_3/\gamma_2$ -HEK
GABA_A $\alpha_3/\beta_3/\gamma_2$ -HEK
GABA_A $\alpha_4/\beta_3/\gamma_2$ -HEK
GABA_A $\alpha_5/\beta_3/\gamma_2$ -HEK
GABA_A $\alpha_6/\beta_3/\gamma_2$ -HEK
GlyRA1-HEK
GluR6-HEK

CRO Services are performed on these cell lines by our sister company Eurofins-Discovery in St. Charles, MO

Human origin unless otherwise noted

Benefits of Ready to Assay (RTA) Cells

- RTA cells eliminate the need for continuous cell culture
- Eliminates inconsistencies that can be inherent in cell line cultures due to:
 - Different cell culture technicians
 - Different time windows of stable cell lines in culture
- RTA cells are frozen down at optimal expression
- Every production lot is QC validated

New PrecisiONRTA cells *Coming soon!*

Cardiac Safety Panel:

- hERG-CHO (CYL3038RTA)
- Nav1.5-HEK (CYL3004RTA)
- Cav1.2 (CYL3051RTA)

Provided with a Certificate of Analysis that includes electrophysiology data and pharmacological characterization for each lot produced

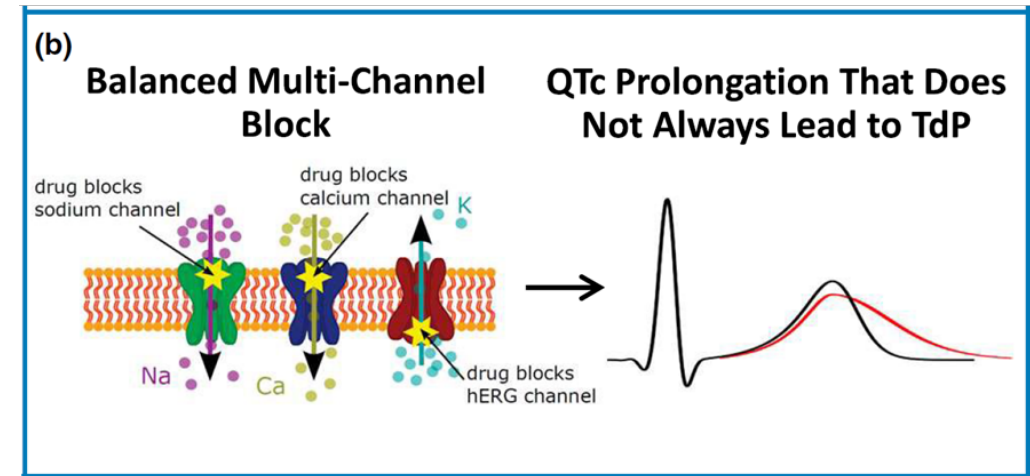
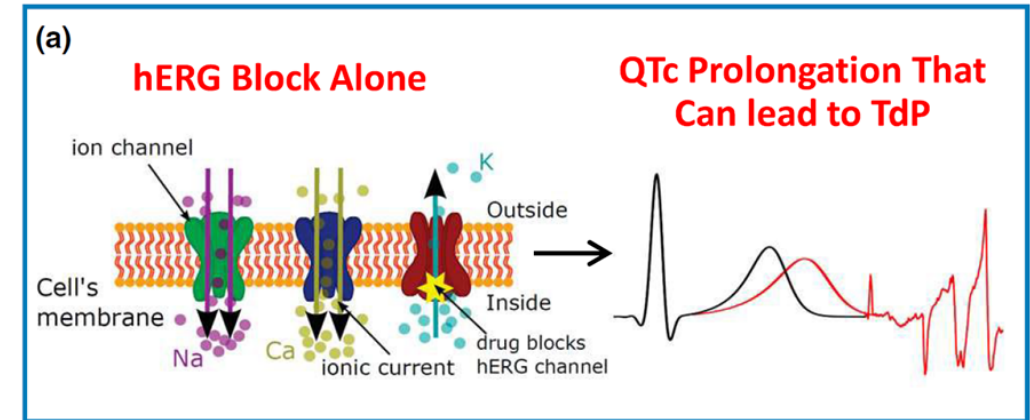
Pain Targets:

- Nav1.8 (CYL3025RTA)
- And other popular therapeutic pain targets

Let us know your favorites!

In support of the new FDA CiPA guidelines, Eurofins DiscoverX® is launching the **Cardiac Safety Panel PrecisiON stable cell lines in a frozen Ready-to-assay (RTA) format**

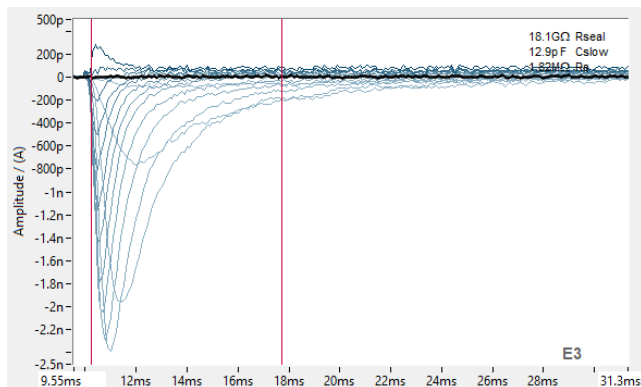
Blockade of Nav1.5 and/or Cav1.2 can offset hERG blockade and render the drug safe.



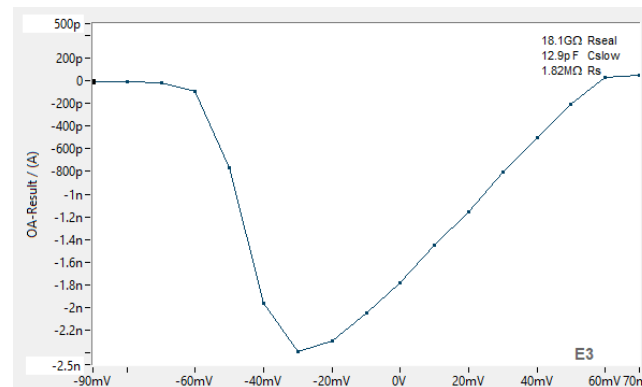
Translational Models and Tools to Reduce Clinical Trials and Improve Regulatory Decision Making for QTc and Proarrhythmia Risk (ICH E14/S7B Updates)

David G. Strauss^{1*}, Wendy W. Wu¹, Zhihua Li¹, John Koerner² and Christine Garnett³

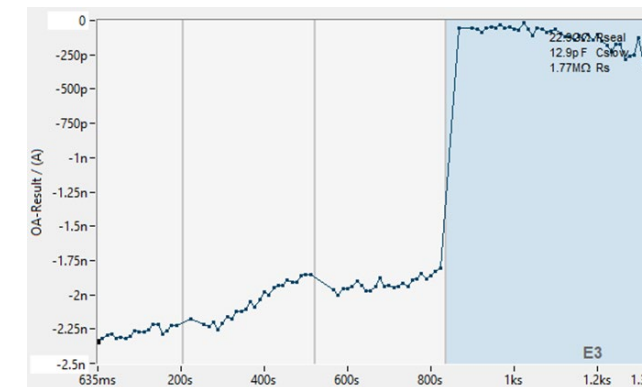
Current Voltage (I/V) Relationship and Tetracaine Pharmacology of PrecisiON™ Nav1.5 Ready-To-Assay Cells



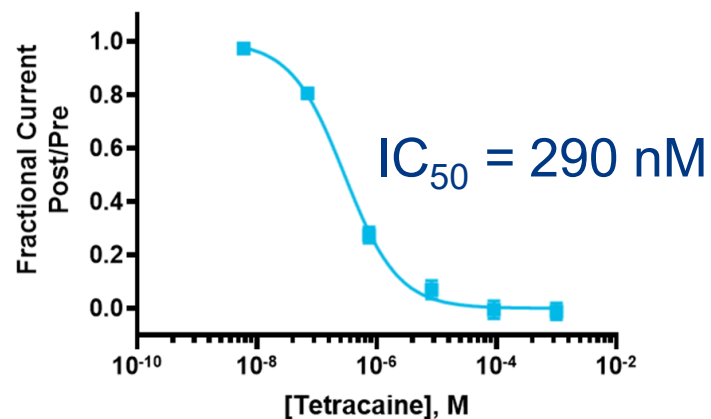
Raw Ionic Currents



I/V Plot



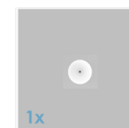
Peak Current over time (I-t plot)



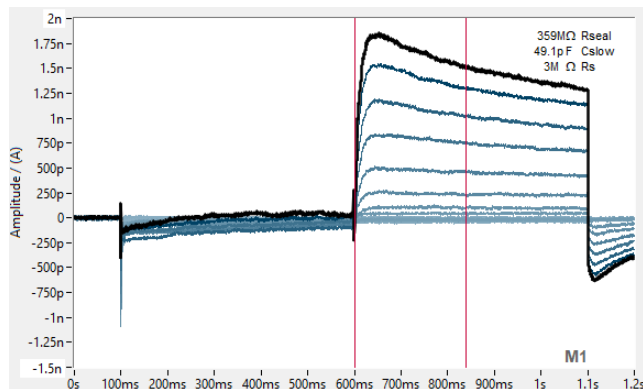
Nav1.5-HEK		
Part #	Channel	Mean Success
CYL3004	Nav1.5	74.0% (n = 2 runs)
CYL3004-RTA	Nav1.5-RTA	71.7% (n = 2 runs)

- Stable RTA currents are obtained that are sensitive to tetracaine
- Success rates are similar stable cell line

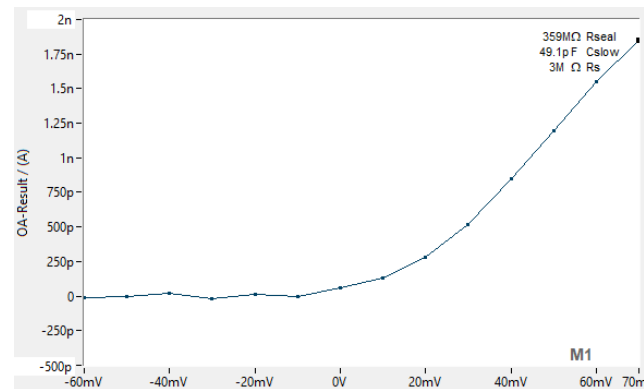
1-Hole/well Recording Plate



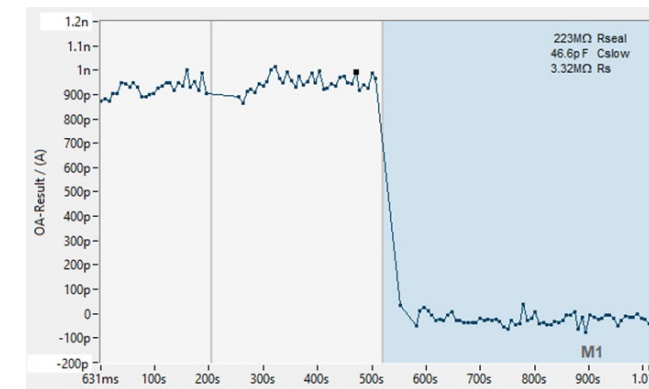
Current Voltage (I/V) Relationship and Cisapride Pharmacology of hERG-CHO Ready-To-Assay Cells



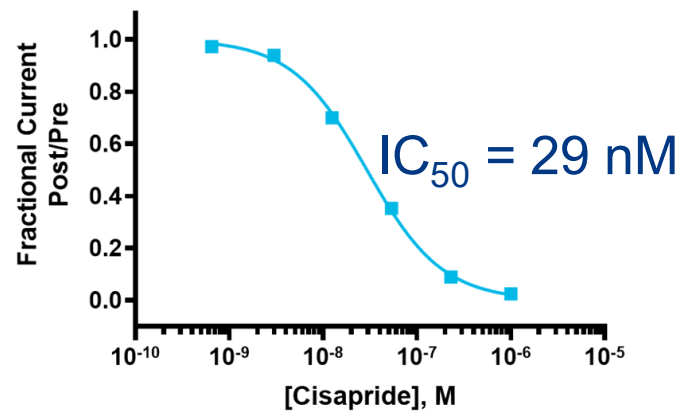
Raw Ionic Currents



I/V Plot



Peak Current over time (I-t plot)



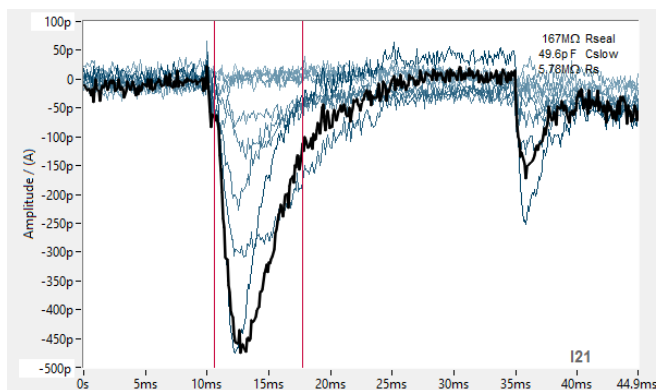
hERG-CHO		
Part #	Channel	Mean Success
CYL3038	hERG Stables	87.5% (n = 2 runs)
CYL3038-RTA	hERG RTA's	83.0% (n = 2 runs)

- Stable RTA currents are obtained that are sensitive to cisapride
- Success rates are similar stable cell line

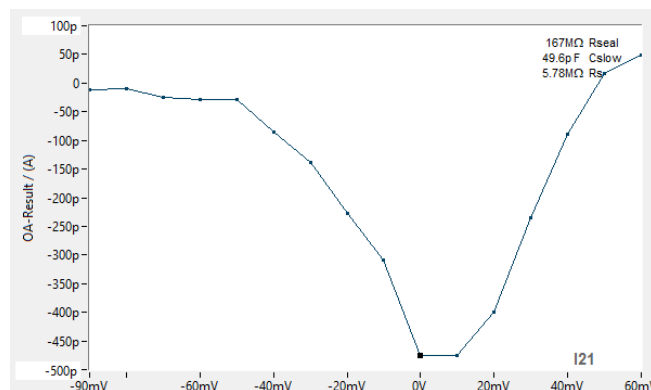
4-Holes/well Population Patch Plate



PrecisION™ Nav1.8 and Cav1.2 Ready-To-Assay Cells

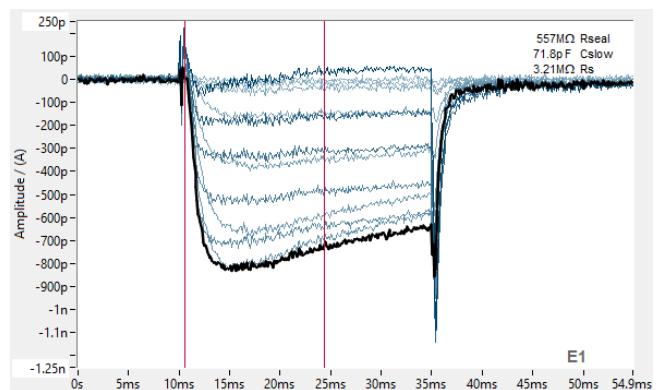


Raw Ionic Currents

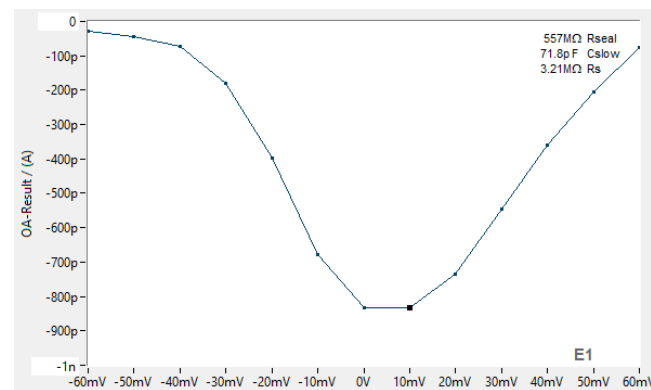


I/V Plot

Nav1.8-HEK		
Part #	Channel	Mean Success
CYL3025-RTA	Nav1.8-RTA	73.2% (n = 2 lots)



Raw Ionic Currents



I/V Plot

Cav1.2-HEK		
Part #	Channel	Mean Success
CYL3051-RTA	Cav1.2-RTA	67.9% (n = 2 lots)

Stable RTA currents are obtained and the success rates are similar stable cell line

- Analgesics designed to inhibit the activity of the Nav1.8 channel in the pain pathway are promising non-addictive alternatives to opioid analgesics
- Latigo has presented compelling evidence that LTGO-33 is a novel Nav1.8 inhibitor
- Eurofins DiscoverX[®] is in the process of converting selected stable cell lines into ready-to-assay format include cardiac safety and therapeutic pain targets



Learn more at discoverx.com/target-class/ion-channel/