



THE ION CHANNEL EXPERT

# Automated patch clamping for screening and profiling in drug discovery and risk assessment

*Emir Duzic*

Ion Channel Retreat, Vancouver  
June 2013

- **ChanTest in Brief**
- **IonWorks Barracuda (IWB) Platform**
  - **Advantages over other platforms**
  - **Assay validation examples:**
    - Nav1.7 validation for screening
    - GABA<sub>A</sub> validation for profiling
    - Cardiac Channel Panel for risk assessment
- **Summary**

# ChanTest at a Glance



- Founded 1998; introduced hERG patch clamp safety assay
- Invented hERG- and Chan-Lite trafficking inhibition assays
- **Highly Trained and Motivated Personnel**
- **~300,000 compounds tested for 350+ clients**
- Successful audits by FDA & 200+ sponsors
- RCA with FDA to improve cardiotox predictivity
- Industry's most complete Ion Channel portfolio
- Recombinant cell lines, SC-derived human cells, primary cells
- Extensive GPCR targets, strong Transporter program
- Named "**most trusted fee-for-service provider**" 2006 – 2012  
(*HTStec Ion Channel Trends Survey*)

# Client Interaction Models



THE ION CHANNEL EXPERT

www.chantest.com



# Discovery Services



THE ION CHANNEL EXPERT

www.chantest.com



<i>Assay Development</i>	<i>Screening</i>	<i>Lead Optimization</i>
<p>Cell lines development &amp; Characterization</p> <p>Custom Assay Development for Primary and Secondary Screening</p>	<p>Assay Ready Cryo-preserved Cells</p> <p>Primary &amp; Secondary Screening Services</p> <p>Hit Triage and Hit to Lead Chemistry</p>	<p><i>In vitro</i> Potency and Efficacy</p> <ul style="list-style-type: none"><li>- Mechanism of Action</li><li>- Selectivity</li><li>- GPCR Target Panels</li></ul> <p>Safety Testing</p> <ul style="list-style-type: none"><li>- hERG Screening Service</li><li>- Disease Channel Panels</li></ul> <p>Lead Optimization Chemistry</p>

**Choose single projects or an entire program**

# Acknowledgements

## ScreenPatch Team

Yuri Kuryshev

Hung Lee

Mike Fraifogl

Cory Stebal

## Channel Pharmacology

Glenn E. Kirsch

## CEO/CSO

Arthur “Buzz” Brown

## Compound Management

Peter Hawryluk

# IonWorks Barracuda Platform

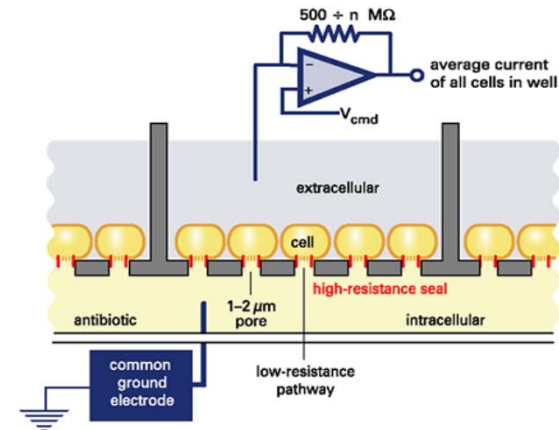


THE ION CHANNEL EXPERT

www.chantest.com

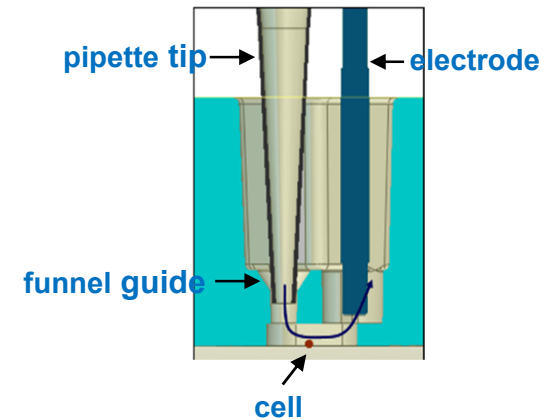


## Planar Patch Method



Cells delivered to electrode

## Flow-through design



- Population Patch Clamp (PPC) and single hole (SH) recording modes (384 well format)
- Controls single cell membrane potential and measures ionic currents in single-cell or cell population ( $\leq 64$  cells/well)
- 384-channel pipettor, integrated 384-channel electronic head
- Features continuous voltage-clamp current measurement with rapid solution addition for fast-desensitizing, ligand-gated channels

# Advantages of IWB Platform



THE ION CHANNEL EXPERT

www.chantest.com

Characteristic	E-Phys Platform		
	IWB	IWQ	QPatch
Continuous voltage clamp	Yes	No	Yes
Seal Resistance	> 300 MΩ	< 50 MΩ	Gigaseal
Parallel recording for all 384 wells	Yes	No	N/A
Possibility of cross-contamination	No	Yes	Yes
Multiple voltage protocols/experiment	Yes	No	Yes
Automated prep of compound plates	Yes <i>(standard)</i>	Yes <i>(standard)</i>	Yes <i>(optional)</i>
Screening capabilities (“Targeted HTS”)	Yes	Yes	No
Manual prep of compound plates	Yes <i>(optional)</i>	Yes <i>(optional)</i>	Yes <i>(standard)</i>
Multiple TA additions	Yes <i>(optional)</i>	No	Yes <i>(standard)</i>
Cost/data point	Moderate	Moderate	High



# Assay Development Process



THE ION CHANNEL EXPERT

www.chantest.com

Cell Line  
Development  
& Validation

*~3 months*

**Expression confirmation:**

- Western blot
- RT-PCR
- Functional cell-based assay

**Cell line optimization:**

- Induction conditions
- Monitor functional expression over 10-12 passages

**Functional Testing:**

- Basic stimulus/response detection
- Test basic functional characteristics
- Reference compounds pharmacology

Assay Development, Optimization  
& Validation

*3 weeks*

- Optimize recording conditions, voltage protocols, intra- & extra-cellular solution (to slow run-down or increase current)
- Adjust stimulus parameters (e.g., ligand addition, voltage steps)
- Stimulus/response testing
- Optimize Cell density
- Optimize SW
- Optimize ZPrime
- DMSO Tolerance
- Signal Stability
- Plate Uniformity
- Intra- and Inter-day variability
- Reference compound pharmacology
- Minimum Significance Ratio (MSR) study

Assay  
Automation

*2 weeks*

- Optimize liquid and plate handling
- Mock Run
- 3 Days Uniformity
- 3 Days MSR tests
- Pilot Screen (10 plates; 3200 cpds)

# Assay Acceptance Criteria - IWB

## Well-level

Parameter	Value
Seal Resistance ( $R_{\text{seal}}$ )	> 100 M $\Omega$ *
Current Amplitude	> 0.2 nA
$R_{\text{seal}}$ Stability	< 50% decrease
Current Stability	< 30% decrease
Voltage Clamp Quality	Visual control

\* $R_{\text{seal}}$  >300 M $\Omega$  for  $\text{Na}_v$ ,  $\text{K}_v$  and  $\text{Ca}_v$  channels

# Assay Acceptance Criteria -IWB

## Plate-level

Parameter	Value
Z' Factor	$\geq 0.4$
CV for MAX Control	$\leq 20\%$
CV for MID Control	$\leq 20\%$
Signal Window (SW)	$\geq 2$
IC <sub>50</sub> for reference antagonists	$\leq 0.5 \log^*$
Voltage Offset (mV)	$< 10 \text{ mV}^{**}$
Current Stability	$< 30\% \text{ decrease}$
Success rate (percent valid wells)	$\geq 90\%$

\* *from historical mean*

\*\* *measured as I-V curve shift*

# Assays available on IWB Platform



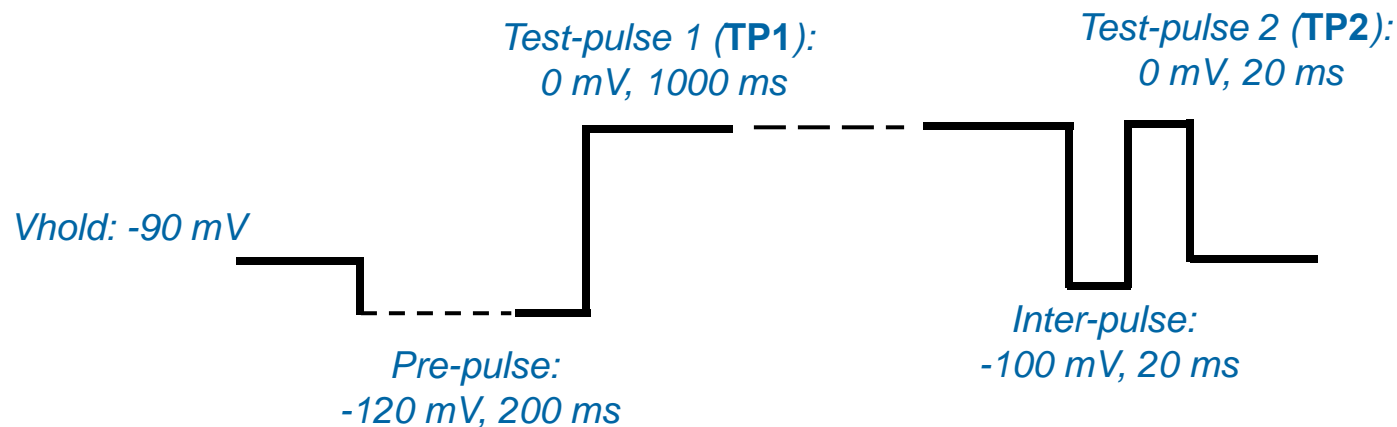
THE ION CHANNEL EXPERT

www.chantest.com

	Ligand Gated Channels	Voltage Gated Ca <sub>v</sub> Channels	Voltage Gated K <sub>v</sub> channels	Voltage Gated Na <sub>v</sub> channels
1	5-HT3A	<b>Cav1.2</b> (β2/cardiac L-type)	hERG	Nav1.1
2	ASIC-1 (ASIC-1a)	Cav2.1 (β4 P/Q type)	Kir2.1	Nav1.2
3	GABA <sub>A</sub> (α <sub>1</sub> β <sub>3</sub> γ <sub>2</sub> )	Cav2.2 (β3/α <sub>2</sub> δ N-type)	Kv1.1	Nav1.3
4	<b>GABA<sub>A</sub></b> (α <sub>2</sub> β <sub>3</sub> γ <sub>2</sub> )	Cav3.2 (T-type)	Kv1.3	Nav1.4
5	GABA <sub>A</sub> (α <sub>3</sub> β <sub>3</sub> γ <sub>2</sub> )		Kv1.4	Nav1.5
6	GABA <sub>A</sub> (α <sub>4</sub> β <sub>3</sub> γ <sub>2</sub> )		Kv1.5	Nav1.6
7	GABA <sub>A</sub> (α <sub>5</sub> β <sub>3</sub> γ <sub>2</sub> )		Kv2.1	<b>Nav1.7</b>
8	HCN2		Kv3.2	Nav1.8
9			Kv3.4	
10			Kv4.3/KChiP2.2	
11			KvLQT1/minK	

# Nav1.7 Assay Protocol on IWB

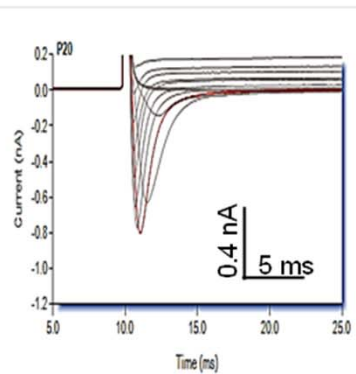
- Cells:
  - CHO cells stably transfected with SCN9A ion channel cDNA (*ChanTest cell line Cat.# CT6003*)
  - Plated 40 - 56 hours before experiments
- Voltage Protocol:



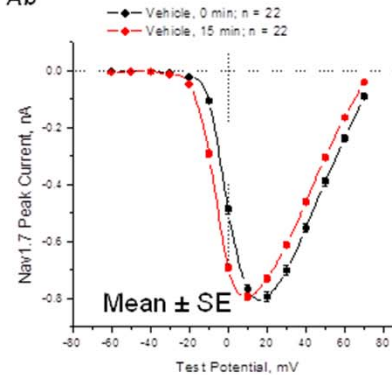
# Nav1.7 Biophysical Characteristics: IWB (PPC) vs. MPC

## A. IonWorks Barracuda

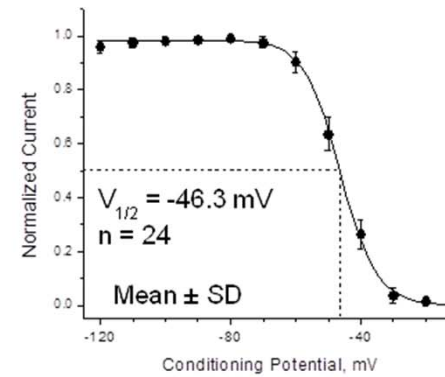
Aa



Ab

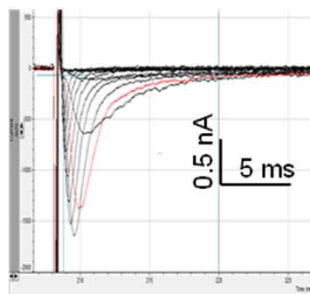


Ac

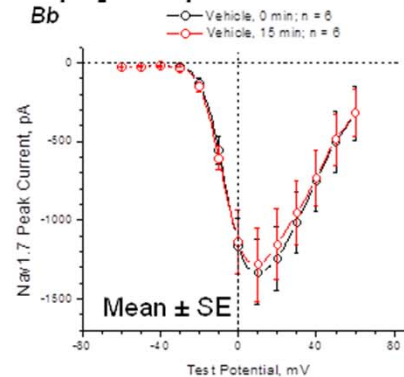


## B. Manual Patch Clamp [Axopatch 200B]

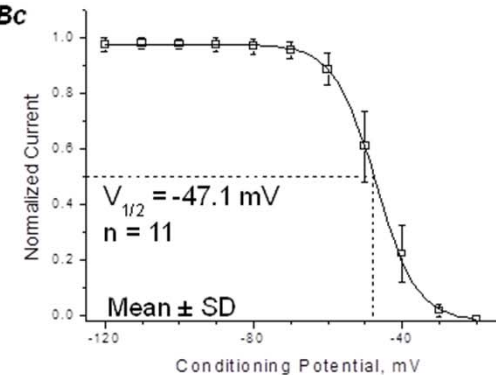
Ba



Bb



Bc



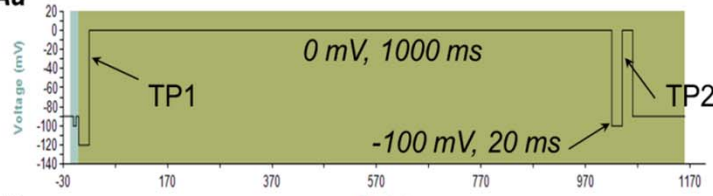
### Voltage protocol:

- holding potential -90 mV.
- 1-second conditioning potential from -120 mV to 50 mV
- 10 mV increment followed by a 20-ms test potential to 0 mV

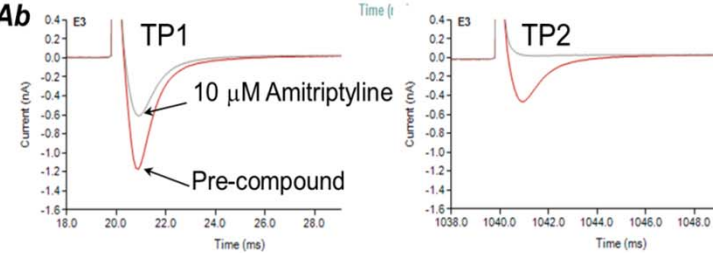
# Inhibition with Amitriptyline: IWB (PPC) vs. MPC

## A. IWB

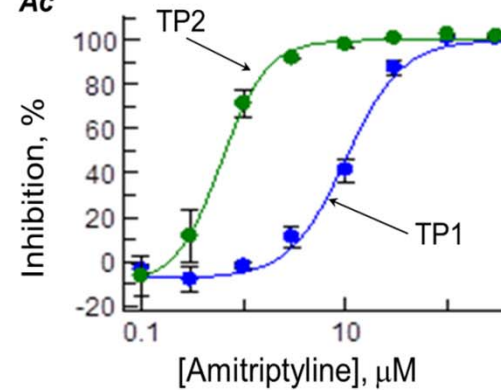
Aa



Ab



Ac



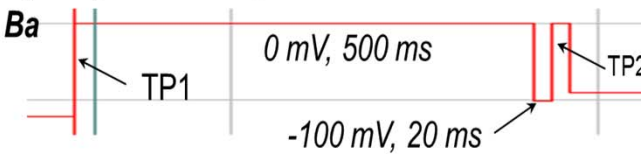
$IC_{50}$  ( $\mu$ M)

TP1 10.01

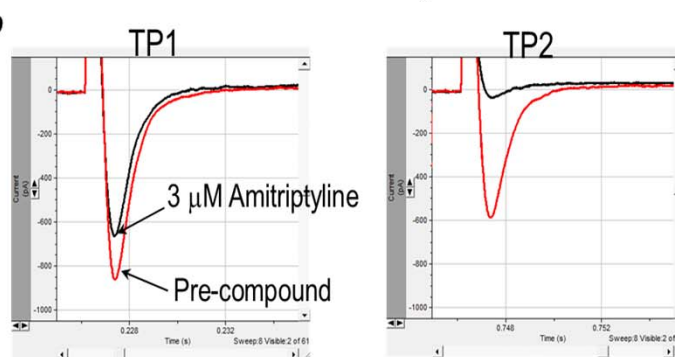
TP2 0.59

## B. MPC [Axopatch 200B]

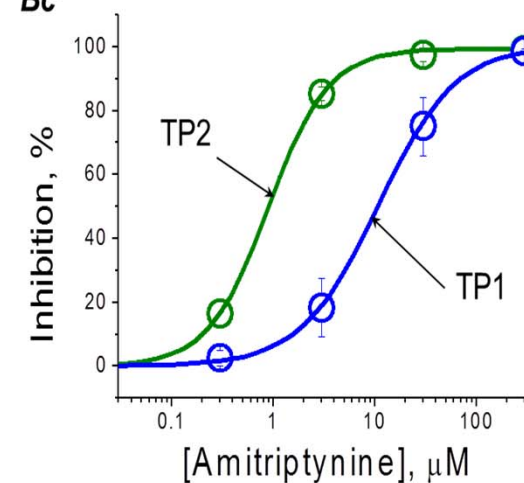
Ba



Bb



Bc



$IC_{50}$  ( $\mu$ M)

TP1 10.80

TP2 0.89

# Reference Compounds Potency: IWB (PPC) vs. MPC

Compound	IWB (IC <sub>50</sub> , μM)*		MPC (IC <sub>50</sub> , μM)**	
	TP1 (tonic block)	TP2 (phasic block)	TP1 (tonic block)	TP2 (phasic block)
Amitriptyline	10.01	0.59	10.80	0.89
Mexiletine	121.94	30.36	116.25	31.46
Flecainide	40.92	16.26	25.07	16.19
Lamotrigine	>300	19.82	631.24	60.15
Carbamazepine	245.70	110.78	659.10	106.30

\*Mean value [8-point CRC (4 replicas/conc.)3 independent exp. (1 exp./day)]

\*\*Mean value [1-2 days/compound; 4-5 CRC (5-8 replicas/conc.)]



# Summary of Uniformity Study



THE ION CHANNEL EXPERT

www.chantest.com

Day	Plate	Mean	SD	%CV	Z'	SW	Success Rate (%)
DAY 1	Max 1	111.68	9.92	8.88	0.651	6.95	88.0
	Max 2	99.79	8.56	8.58			89.3
	Mid 1	57.27	7.22	12.61			83.6
	Mid 2	62.46	7.32	11.71			87.0
	Min 1	6.51	2.20	33.74			91.9
	Min 2	7.04	2.31	32.75			91.1
DAY 2	Max 1	91.22	9.94	10.90	0.566	4.85	88.8
	Max 2	95.36	10.05	10.54			91.9
	Mid 1	47.75	7.08	14.82			89.8
	Mid 2	57.45	8.46	15.04			93.3
	Min 1	5.85	2.38	40.80			89.3
	Min 2	9.23	2.48	26.85			92.7
DAY 3	Max 1	94.89	11.47	12.08	0.586	5.14	82.0
	Max 2	100.38	9.29	9.25			83.3
	Mid 1	49.51	8.57	17.20			85.7
	Mid 2	61.14	8.71	14.25			84.9
	Min 1	6.56	2.85	43.46			79.9
	Min 2	5.70	1.23	21.56			85.9

Max signal (Vehicle); Mid Signal (150  $\mu$ M Mexiletine); Min Signal (1000  $\mu$ M Mexiletine)

# Mexiletine Uniformity at TP1 and TP2



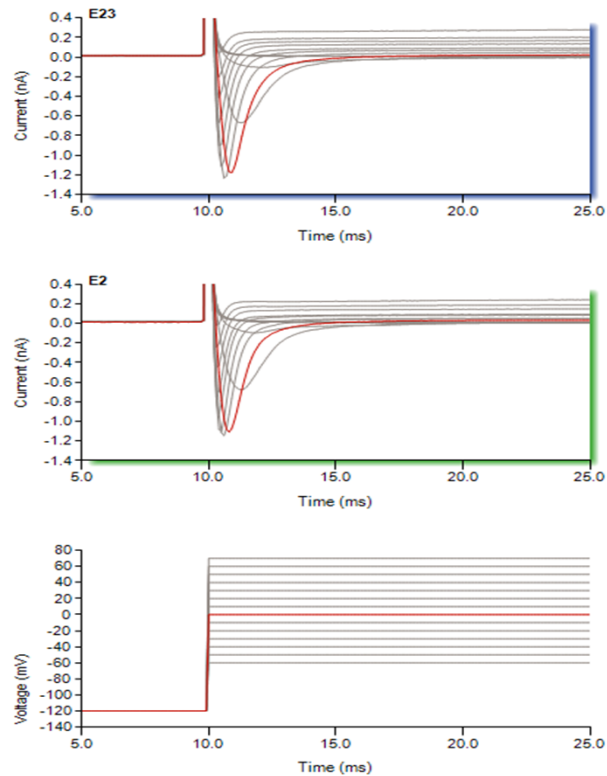
THE ION CHANNEL EXPERT

www.chantest.com

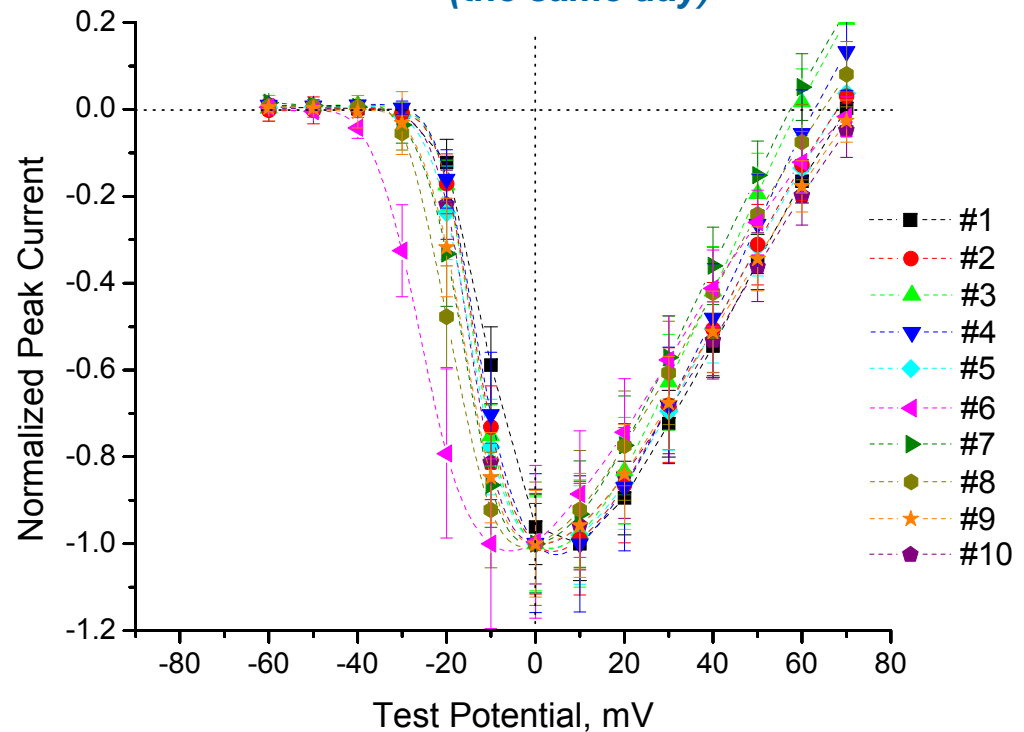
Experiment ID	Z'		Mexiletine IC50 (μM)		Success Rate (%)
	T1	T2	T1	T2	
Day 1 #1	0.63	0.63	98.5	23.1	83.9
Day 1 #2	0.71	0.82	163.2	28.2	90.1
Day 1 #3	0.67	0.68	73.2	9.5	88.8
Day 1 #4	0.67	0.71	128.0	19.7	90.1
Day 1 #5	0.67	0.51	67.6	16.1	79.7
Day 1 #6	0.56	0.43	120.3	13.3	80.7
Day 1 #7	0.71	0.68	123.8	16.6	89.6
Day 1 #8	0.51	0.58	136.0	20.3	96.6
Day 1 #9	0.67	0.67	159.1	36.2	94.8
Day 1 #10	0.66	0.68	246.6	37.6	96.1
Day 2 #1	0.81	0.78	105.0	43.8	96.4
Day 3 #1	0.76	0.79	162.4	24.1	95.3
<b>Mean</b>	<b>0.67</b>	<b>0.66</b>	<b>132.0</b>	<b>24.0</b>	<b>90.2</b>
<b>SD</b>	<b>0.08</b>	<b>0.11</b>	<b>48.1</b>	<b>10.5</b>	<b>6.1</b>
<b>Range</b>	<b>0.51 - 0.81</b>	<b>0.51 - 0.82</b>	<b>67.6 - 246.6</b>	<b>9.5 - 43.8</b>	<b>79.7 - 96.6</b>

# Daily Stability of Voltage Potentials

## A. Representative Nav1.7 current traces



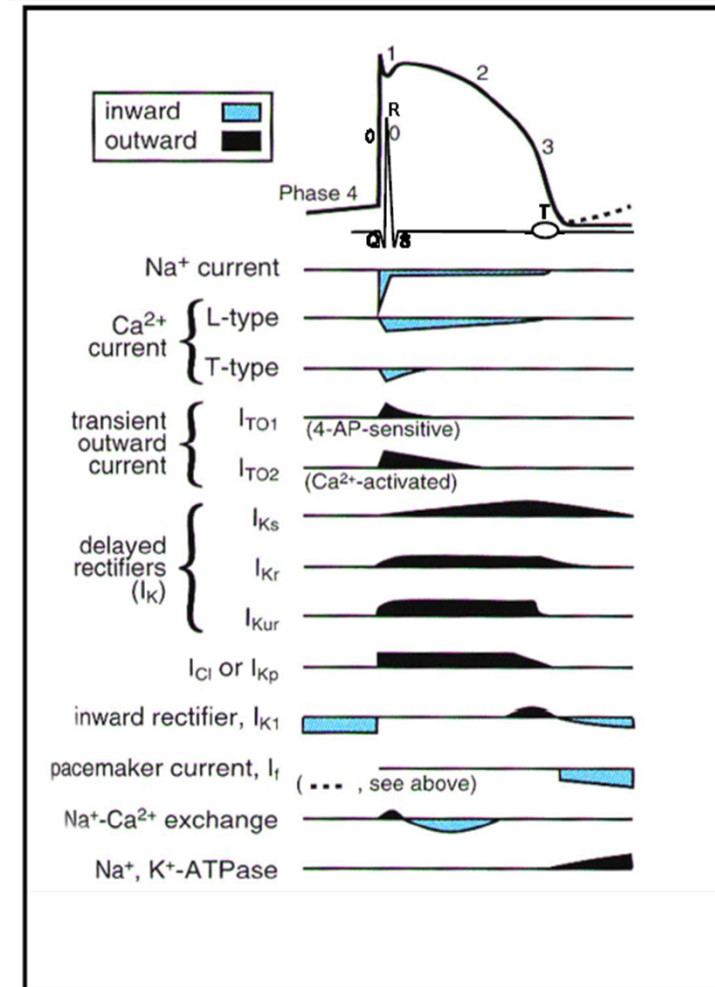
## B. I-V curves for 10 independent experiments (the same day)



Data presented as Mean  $\pm$  SD ( $n = 26 - 32$ ).

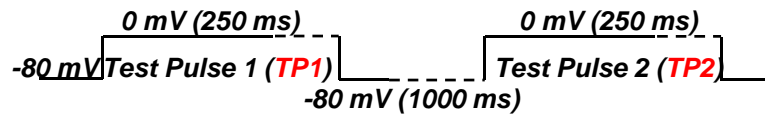
# It is Not Just About hERG

Cardiac Channel Panel™	Action Potential Phase
Nav1.5 ( $I_{Na}$ )	0,2
Cav 1.2/ $\beta$ 2, $\alpha$ 2 $\delta$ , (L-type)	2
Cav 3.2 (T-type)	1
Kv4.3 ( $I_{TO1}$ )	1
KvLQT1/minK ( $I_{KS}$ )	2-3
hERG ( $I_{Kr}$ )	2-3
Kv1.5 ( $I_{Kur}$ )	2-3
Kir2.1 ( $I_{K1}$ )	4
HCN2 (pacemaker, $I_f$ )	4
HCN4 (pacemaker, $I_f$ )	4
Kir3.1/3.4 ( $I_{K,Ach}$ )	4
Kir6.2/SUR2A ( $I_{K,Ach}$ )	4
NCX1 (Na-Ca exchange)	2

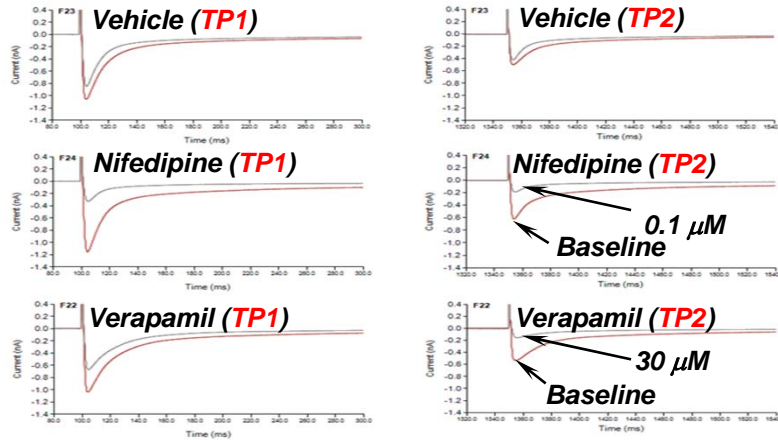


# Cardiac Channel Panel: Cav1.2/ $\beta$ 2/ $\alpha$ 2 $\delta$ 1 Protocol on IWB: Uniformity of Nifedipine and Verapamil Inhibition

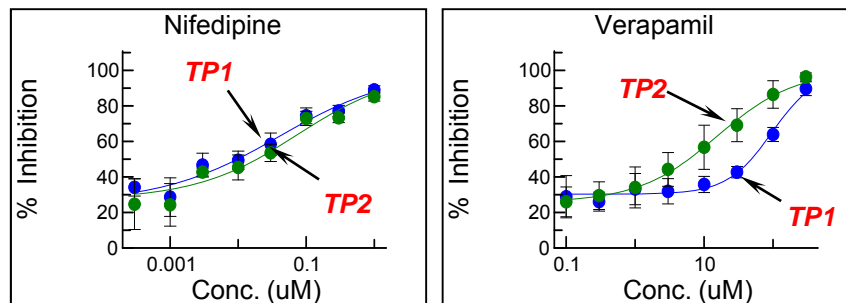
## A. Voltage Protocol



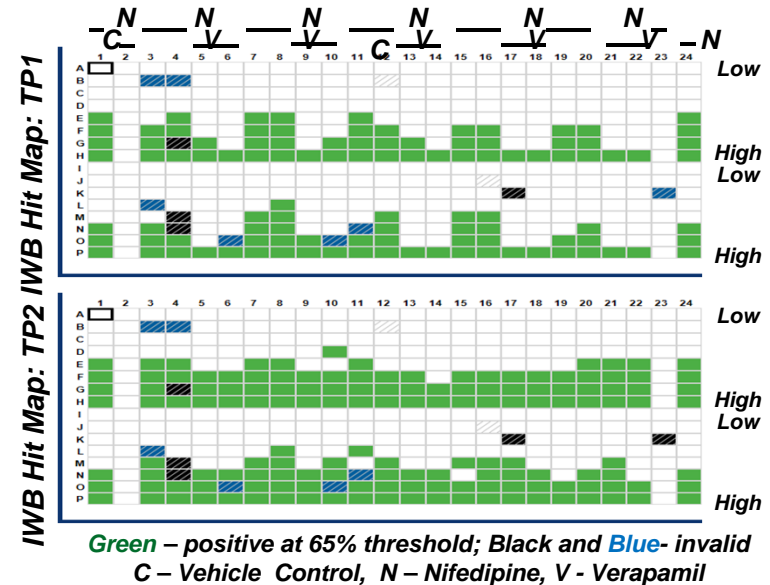
## B. Representative Current Traces



## C. Representative Dose-Response Curves



## D. Hit Map



## E. Intra-plate Uniformity

### Nifedipine Block

Column	IC <sub>50</sub> ( $\mu$ M)	
	TP1	TP2
3, 4	0.044	0.074
7, 8	0.028	0.035
11, 12	0.045	0.067
15, 16	0.035	0.043
19, 20	0.036	0.058
Mean	0.038	0.055
SD	0.007	0.016
Range	0.028-0.045	0.035-0.074

### Verapamil Block

Column	IC <sub>50</sub> ( $\mu$ M)	
	TP1	TP2
5, 6	96.8	15.4
9, 10	81.7	13.3
13, 14	115.4	14.4
17, 18	81.5	14.0
21, 22	118.1	19.1
Mean	98.7	15.2
SD	17.6	2.3
Range	81.5-118.1	13.3-19.1

# Example of Cardiac Panel Report



THE ION CHANNEL EXPERT

www.chantest.com

Program Code	Study Submission ID	IC <sub>50</sub> , μM											
		Cav1.2		Nav1.5		hERG	Kv1.5		Kv4.3/KChIP2.2		LQT1/ minK	Kir2.1	HCN2
		TP1	TP2	TP1	TP2	Peak	Peak	300-ms	Peak	25-ms	2000-ms	500-ms	2000-ms
1	CPD_00845_02_A	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
2	CPD_00854_02_A	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
3	CPD_00872_01_A	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
4	CPD_00905_01_A	>30	24.73	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
5	CPD_00822_03_A	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
6	CPD_00745_03_A	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
7	CPD_00801_02_A	26.03	12.43	>30	>30	>30	23.57	17.05	>30	29.35	>30	>30	>30
8	CPD_00763_02_A	>30	>30	>30	>30	>30	>30	23.66	>30	>30	>30	>30	>30
9	CPD_00449_01_A	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30	>30
10	CPD_00524_01_A	>30	>30	>30	>30	>30	>30	25.16	>30	14.92	>30	>30	>30
PC	Nifedipine	0.050	0.055										
PC	Lidocaine			1035.92	12.92								
PC	E-4031					0.032							
PC	4-AP						1173.74	97.92					
PC	Flecainide								61.13	6.57			
PC	Chromanol										3.14		
PC	Ba <sup>2+</sup>											7.96	
PC	ZT-7288												11.35

# Summary

- Using IonWorks Barracuda (IWB) platform we have developed 31 robust electrophysiology assays for screening and profiling compounds against voltage-gated and ligand-gated ion channels
- All assays were validated using strict acceptance criteria at the well and plate level
- For voltage-gated channels assay protocols has been configured to assess both tonic and use dependent block
- Assays validated on IWB platform can be used at early stage of drug discovery process for rapid cardiovascular safety risk assessment



THE ION CHANNEL EXPERT

***Thank You for your attention!***

***Emir Duzic***

**[eduzic@chantest.com](mailto:eduzic@chantest.com)**

***Tel 216 584-0531***

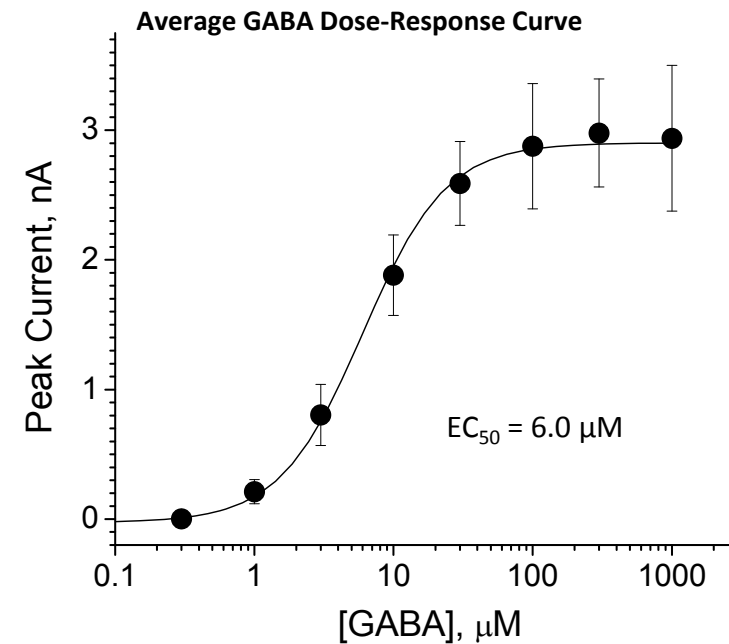
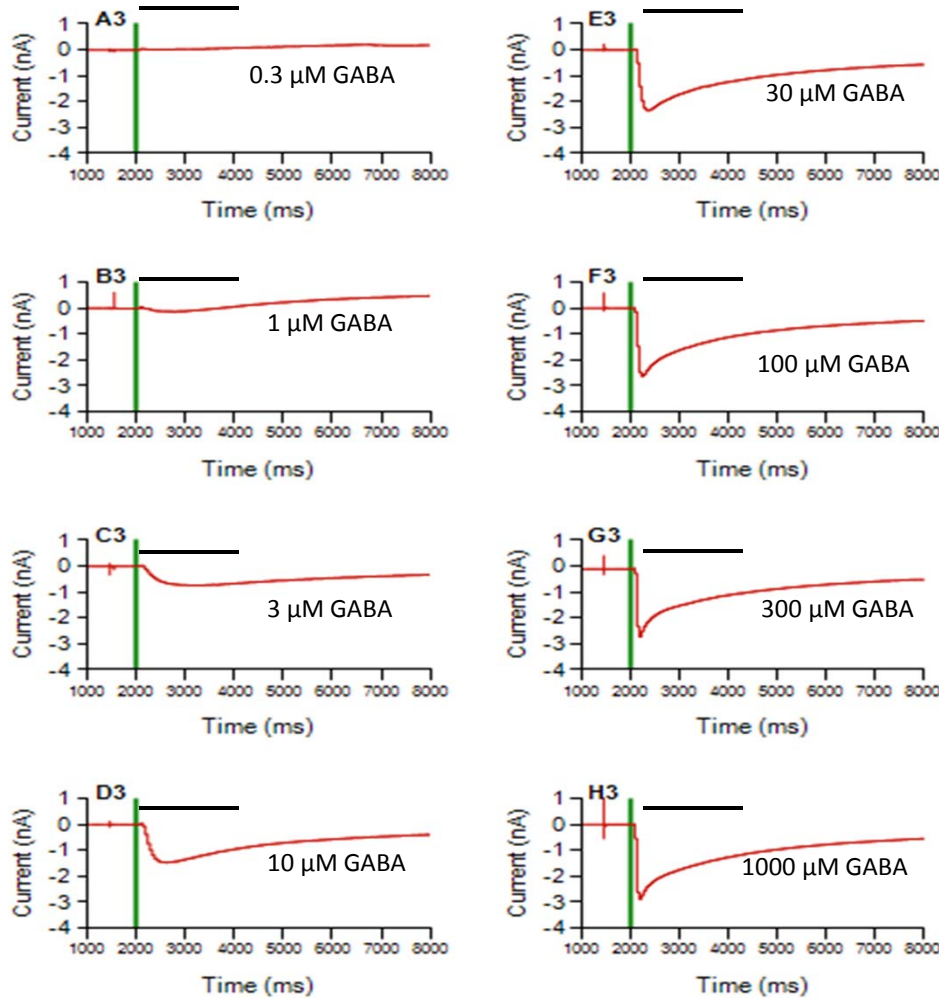




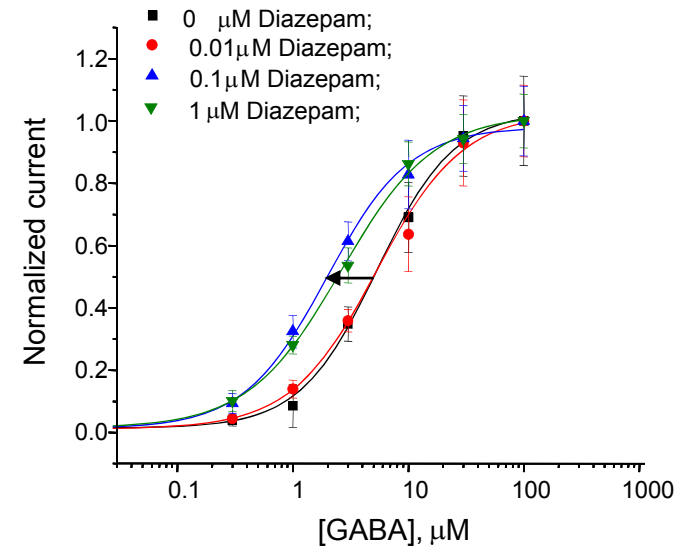
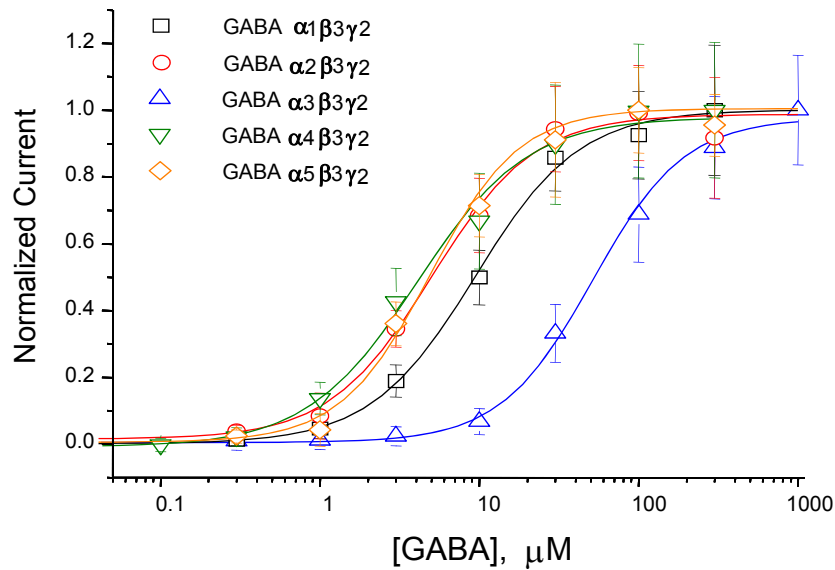
# Backup

# GABA<sub>A</sub> $\alpha_2\beta_3\gamma_2$ -HEK: GABA Concentration-Response

GABA-induced Currents (Wells A3 – H3)



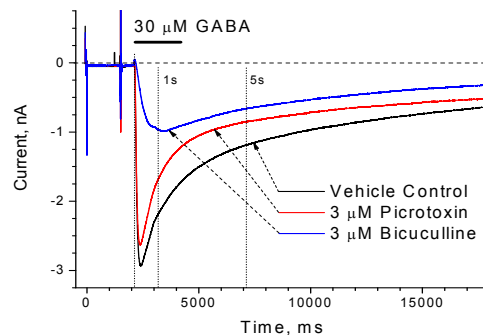
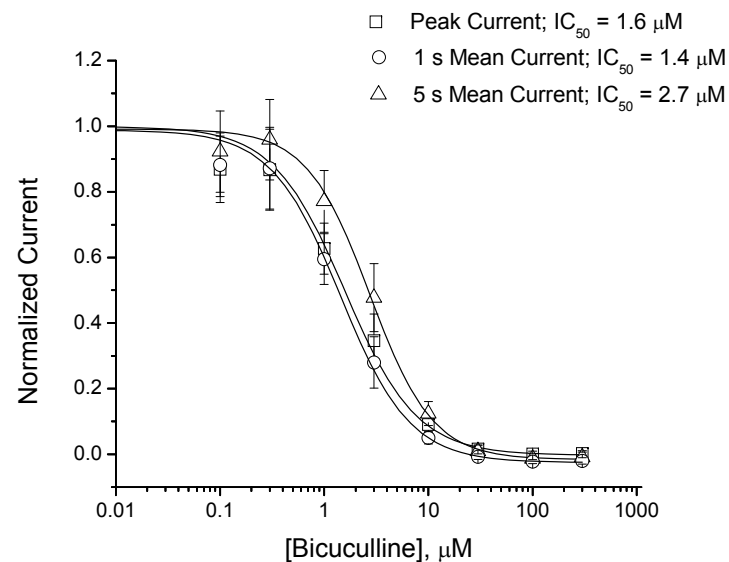
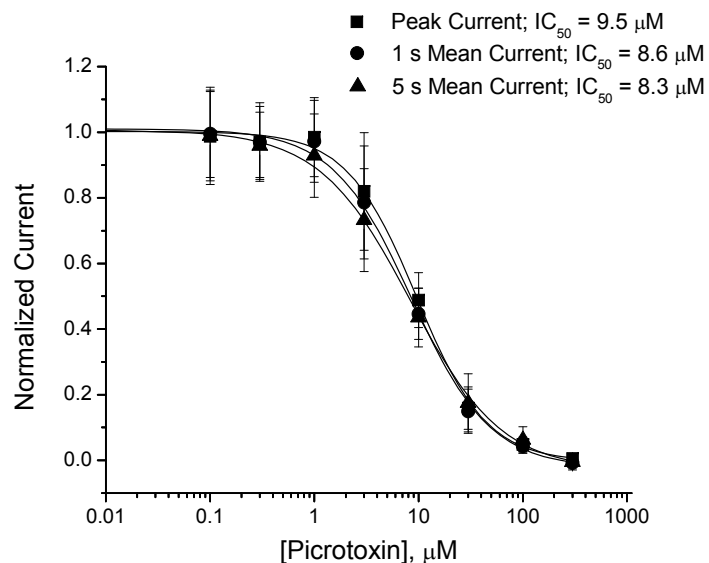
# GABA<sub>A</sub>( $\alpha_x\beta_3\gamma_2$ ) Profiling: Agonist and PAM Modes



GABA <sub>A</sub> Receptor	EC <sub>50</sub> (μM)			
	Mean	SD	N	Publish Values
$\alpha 1/\beta 3/\gamma 2$	12.7	4.2	4	8.3
$\alpha 2/\beta 3/\gamma 2$	5.4	0.7	4	
$\alpha 3/\beta 3/\gamma 2$	47.1	20.3	4	34.7
$\alpha 4/\beta 3/\gamma 2$	4.1	0.3	4	3.9
$\alpha 5/\beta 3/\gamma 2$	4.7	0.8	3	

Diazepam (μM)	GABA EC <sub>50</sub> (μM)
0	5.4
0.01	5.3
0.1	2.0
10	2.6

# GABA<sub>A</sub> ( $\alpha_2\beta_3\gamma_2$ ) Profiling: Antagonist Mode



Superimposition of representative traces of the GABA-induced current