

# QPatch Project Report

Generated 2015-04-16 14:40 by KTS

## Project Summary

Project name: Nav inactivation experiment  
Project ID: 1316p  
Database: qdb on 192.168.20.31, version 5.2.0.30527  
Project notes: Demonstration report 1  
Nav inactivation experiment

Last analysis modification: 2015-04-16 14:26  
Last data modification: 2015-04-16 14:26

## Job Summaries

**Job name:** KTA\_Nav inactivation **Job ID: 812**  
**Assay name:** Nav inactivation  
**Created:** 2013-10-25 10:05 by KTA  
**Executed:** 2013-10-25 15:40 by YSH  
**Screening station:** QPATCH255, SW version 4.1.2 (16449) Falster  
**Status:** FINISHED  
**Pipetting:** Voltage gated **Dose-response job:** Yes  
**QPlate type:** QPlate Large **Compound plate format:** MTP-96 8x(3x4)  
**Intracellular solution:** IC. 5.0.0 **Extracellular solution:** EC 0.0.0  
**Compound list:** tonic and use tetracaine 4DR  
**Experiments completed:** 15 **Experiments started:** 15  
**Used compound plates:** 1 **Bar code(s):** 00000  
**Used QPlates:** 1 **Bar code(s):** 01075130013441  
**Success rates:** Priming: 100.0%, Cell positioning: 93.8%, Gigaseal: 93.8%, Whole-cell: 93.8%  
**Job notes:** NaCl50

## Assay Summaries

**Assay name:** Nav inactivation  
**Cell type:** HEK293 **Ion channel:** Nav1.7  
**Whole-cell protocol:** \*HEK Standard  
**Application protocol:** Nav slow inactivation  
**Voltage protocol:** steady state inactivation

## Experiment Results

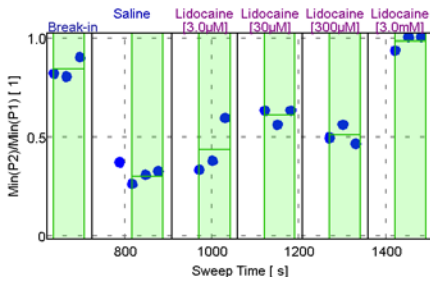
Experiment	XC50(HillFit) [ $\mu$ M]	NoPnt(HillFit)	FitQual(HillFit)	
Exp 812.4.1 (Lidocaine)		3.46m	4	0.06
Exp 812.4.2 (Lidocaine)		50.92m	4	0.03
Exp 812.5.1 (Lidocaine)		3.10m	4	0.01
Exp 812.5.2 (Lidocaine)		260.51 $\mu$	4	0.01
Exp 812.6.1 (Lidocaine)		2.75 $\mu$	4	0.03
Exp 812.1.1 (Tetracaine)		2.77 $\mu$	4	0.03
Exp 812.1.2 (Tetracaine)		1.98 $\mu$	4	0.00
Exp 812.1.3 (Tetracaine)		2.69 $\mu$	4	0.03
Exp 812.2.1 (Tetracaine)		1.73 $\mu$	4	0.00
Exp 812.2.2 (Tetracaine)		3.04 $\mu$	4	0.00
Exp 812.2.3 (Tetracaine)		2.27 $\mu$	4	0.00
Exp 812.3.1 (Tetracaine)		2.62 $\mu$	4	0.00
Exp 812.3.2 (Tetracaine)		3.61 $\mu$	4	0.00
Exp 812.3.3 (Tetracaine)		3.37 $\mu$	4	0.00

## Experiment Group Results

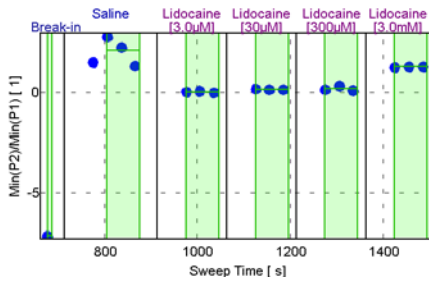
Group Name	XC50(GrpHillFit) [ $\mu$ M]	Avg(XC50(HillFit)) [ $\mu$ M]	NoPnt(GrpHillFit)	FitQual(GrpHillFit)
Group Lidocaine	666.77m	11.55m	20	1.05
Group Tetracaine	2.31 $\mu$	2.68 $\mu$	36	0.51

Xt-plot, input: Sweep Time, Min(P2)/Min(P1)

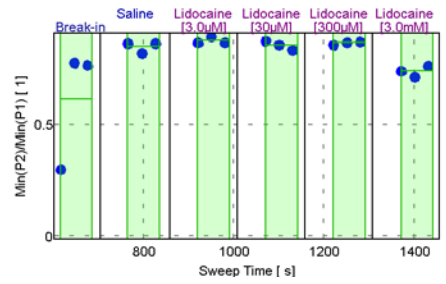
Exp 812.4.1 (Lidocaine)



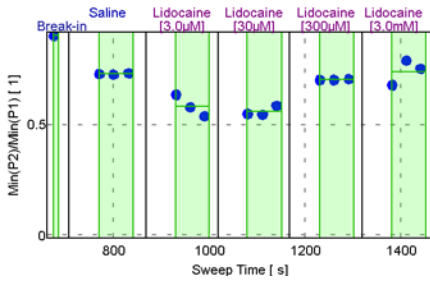
Exp 812.4.2 (Lidocaine)



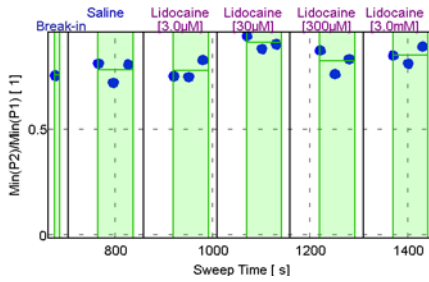
Exp 812.5.1 (Lidocaine)



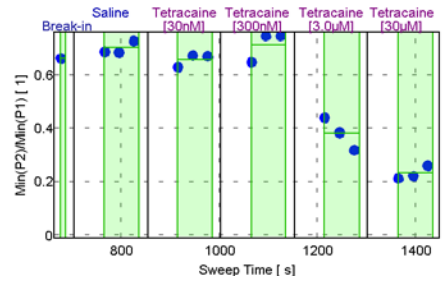
Exp 812.5.2 (Lidocaine)



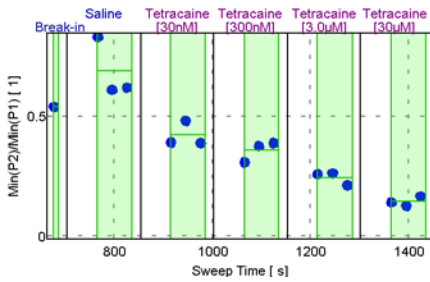
Exp 812.6.1 (Lidocaine)



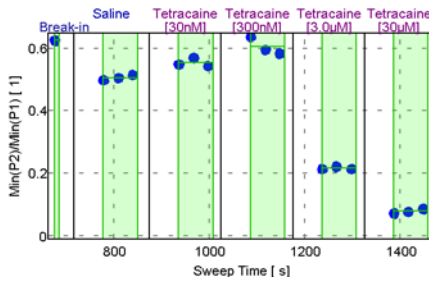
Exp 812.1.1 (Tetracaine)



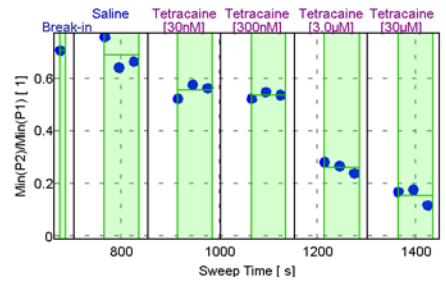
Exp 812.1.2 (Tetracaine)



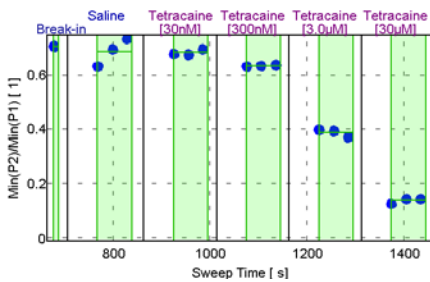
Exp 812.1.3 (Tetracaine)



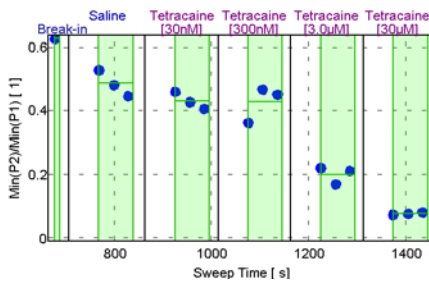
Exp 812.2.1 (Tetracaine)



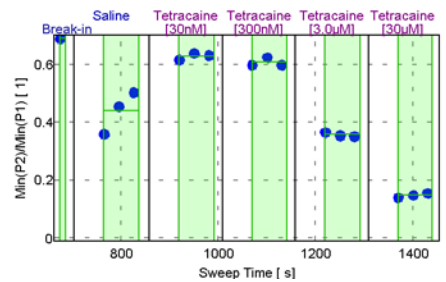
Exp 812.2.2 (Tetracaine)



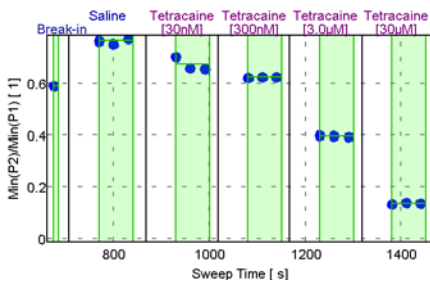
Exp 812.2.3 (Tetracaine)



Exp 812.3.1 (Tetracaine)



Exp 812.3.2 (Tetracaine)



Exp 812.3.3 (Tetracaine)

