

# QPatch Project Report

Generated 2015-04-16 15:23 by KTS

## Project Summary

Project name: KTS\_hERG 5 pt. DR demo  
Project ID: 1319p  
Database: qdb on 192.168.20.31, version 5.2.0.30527  
Project notes: Auto generated from job ID1420 (KTS\_hERG 5 pt. DR demo)  
Based on the template 'KTS\_hERG 5 pt. DR demo 5 uL (id 1420)' which was modified on 2015-04-16 15:16:27

Last analysis modification: 2015-04-16 15:19  
Last data modification: 2015-04-16 15:19

## Job Summaries

**Job name:** KTS\_hERG 5 pt. DR demo **Job ID: 1420**  
Assay name: hERG 5 pt. DR demo 5 uL  
Created: 2015-03-26 12:04 by KTS  
Executed: 2015-03-26 12:10 by KTS  
Screening station: QPATCH262, SW version 5.2.34 (31081) Nyborg  
Status: FINISHED  
Pipetting: Voltage gated Dose-response job: Yes  
QPlate type: QPlate Large Compound plate format: MTP-96 8x(2x6)  
Intracellular solution: IC 0.0.0. Extracellular solution: EC 0.0.0.  
Compound list: hERG validation 5DR 4sticky cmp  
Experiments completed: 15 Experiments started: 20  
Used compound plates: 1 Bar code(s): 00000  
Used QPlates: 2 Bar code(s): 00675140054375,  
00675140054392  
Success rates: Priming: 100.0%, Cell positioning: 93.8%, Gigaseal: 78.1%, Whole-cell: 65.6%  
Job notes:

## Assay Summaries

**Assay name:** hERG 5 pt. DR demo 5 uL  
Cell type: CHO Ion channel: hERG  
Whole-cell protocol: \*CHO 2005-08-17  
Application protocol: hERG 5 pt. DR demo 5 uL  
Voltage protocol: hERG 2 sec activation

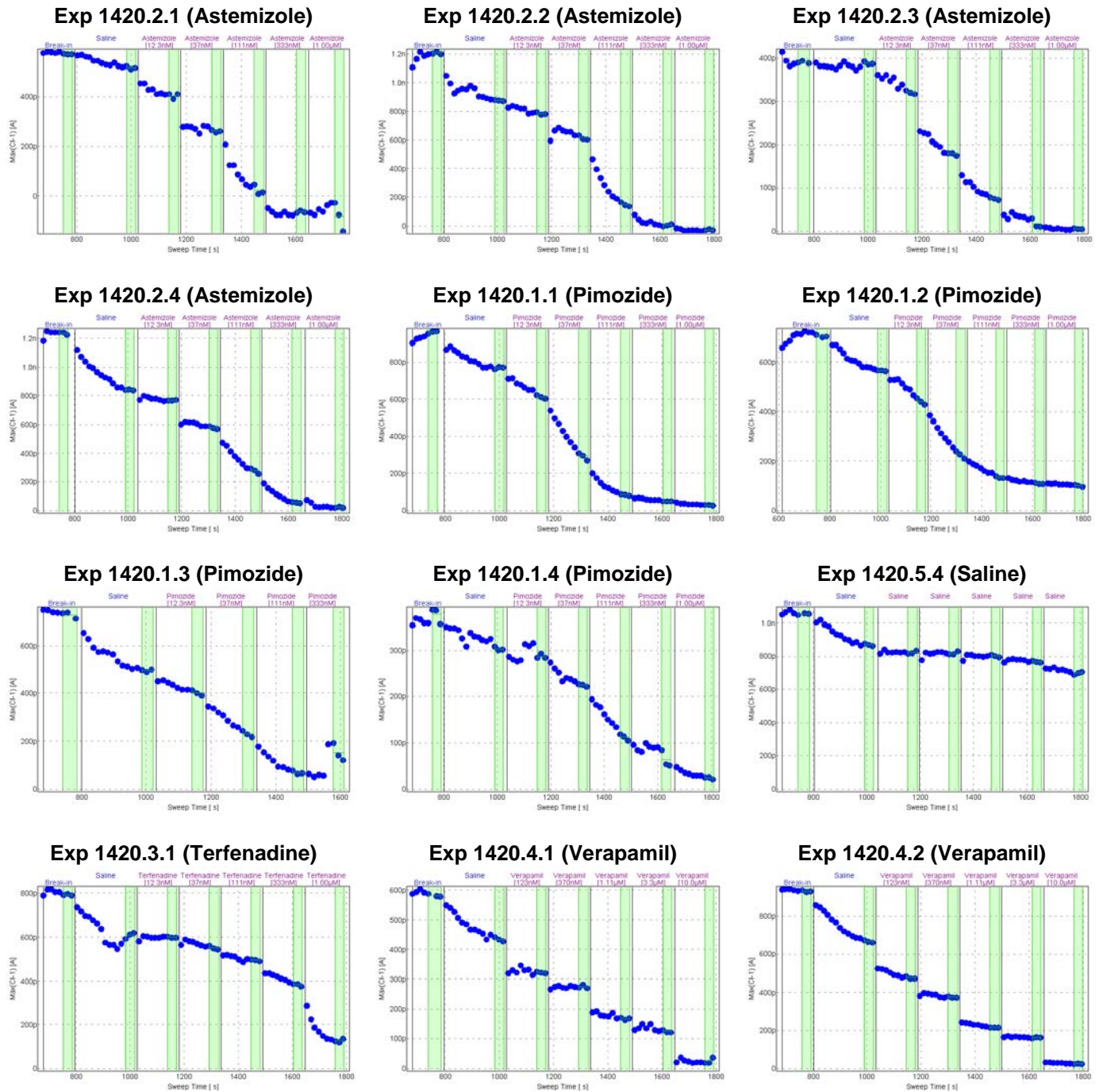
## Experiment Results

Experiment	Sal[Avg(I <sub>t</sub> -plot)] [pA]	FitQual(HillFit)	NoPnt(HillFit)	XC50(HillFit) [μM]	
Exp 1420.2.1 (Astemizole)	516.39p	0.14	5	32.78n	
Exp 1420.2.2 (Astemizole)	877.11p	0.04	5	53.30n	
Exp 1420.2.3 (Astemizole)	388.61p	0.02	5	35.33n	
Exp 1420.2.4 (Astemizole)	842.26p	0.01	5	65.03n	
Exp 1420.1.1 (Pimozide)	770.98p	0.03	5	28.20n	
Exp 1420.1.2 (Pimozide)	566.79p	0.10	5	34.27n	
Exp 1420.1.3 (Pimozide)	496.38p	0.15	4	37.21n	
Exp 1420.1.4 (Pimozide)	304.08p	0.04	5	85.31n	
Exp 1420.5.4 (Saline)	870.47p	Failed	0	Failed	
Exp 1420.3.1 (Terfenadine)	607.53p	0.04	5	421.24n	
Exp 1420.4.1 (Verapamil)	435.26p	0.06	5	658.91n	
Exp 1420.4.2 (Verapamil)	667.46p	0.05	5	469.61n	
Exp 1420.4.3 (Verapamil)	721.32p	0.06	5	582.89n	
Exp 1420.4.4 (Verapamil)	619.12p	0.04	5	412.19n	

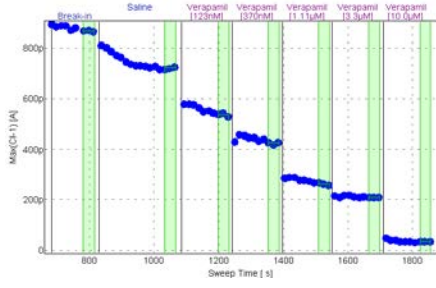
## Experiment Group Results

Group Name	XC50(GrpHillFit) [nM]	FitQual(GrpHillFit)	NoPnt(GrpHillFit)
Group		Failed	Failed
Group Astemizole	44.05n	0.26	20
Group Pimozide	41.85n	0.42	19
Group Saline	Failed	Failed	0
Group Terfenadine	421.24n	0.17	5
Group Verapamil	521.53n	0.20	20

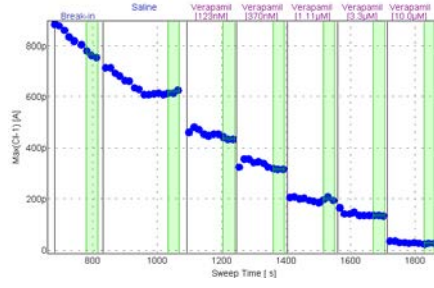
## It-plot, input: Sweep Time, Max(CI-1)



Exp 1420.4.3 (Verapamil)

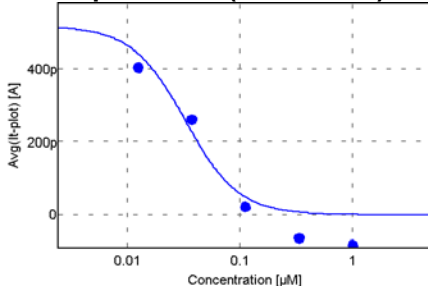


Exp 1420.4.4 (Verapamil)

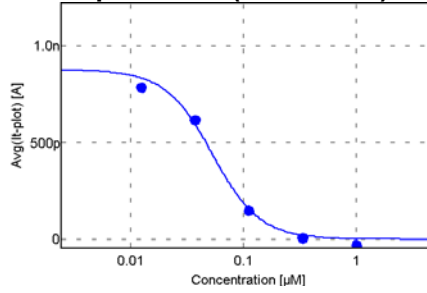


HillFit, input: Concentration, Avg(It-plot)

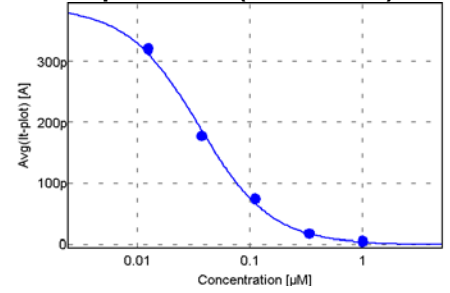
Exp 1420.2.1 (Astemizole)



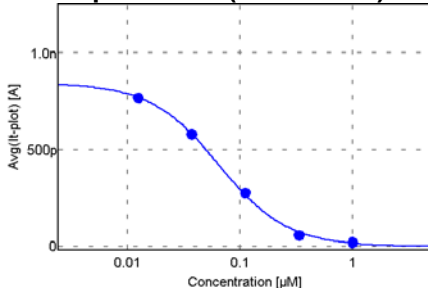
Exp 1420.2.2 (Astemizole)



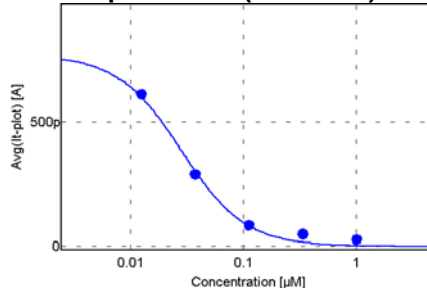
Exp 1420.2.3 (Astemizole)



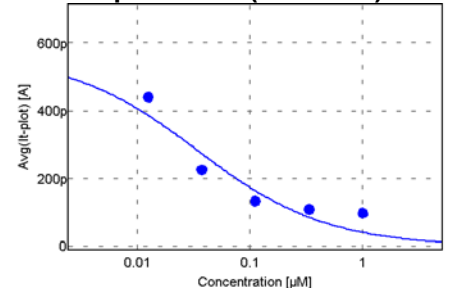
Exp 1420.2.4 (Astemizole)



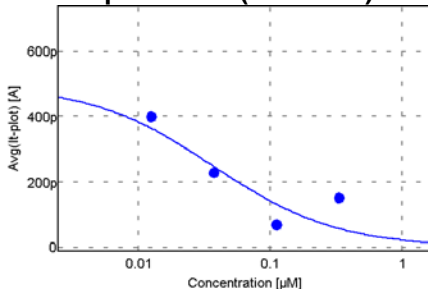
Exp 1420.1.1 (Pimozide)



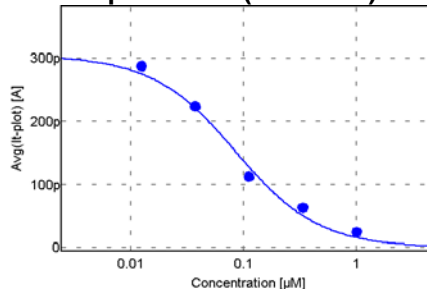
Exp 1420.1.2 (Pimozide)



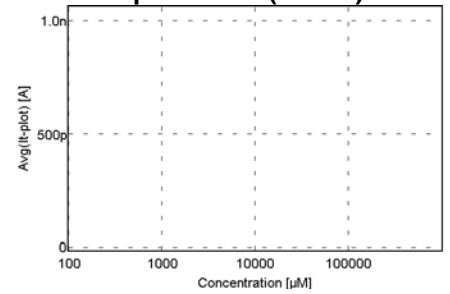
Exp 1420.1.3 (Pimozide)



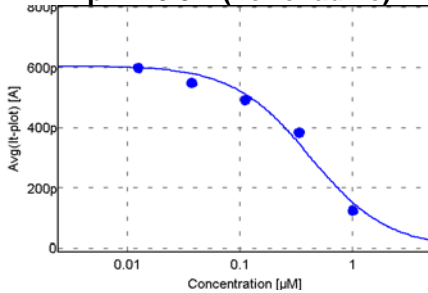
Exp 1420.1.4 (Pimozide)



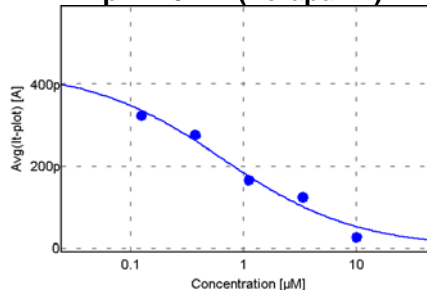
Exp 1420.5.4 (Saline)



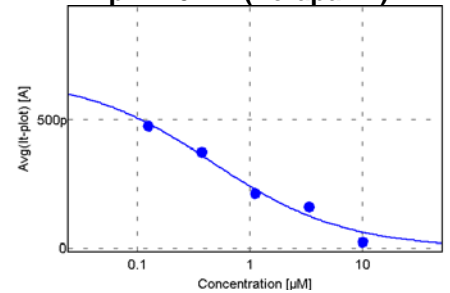
Exp 1420.3.1 (Terfenadine)



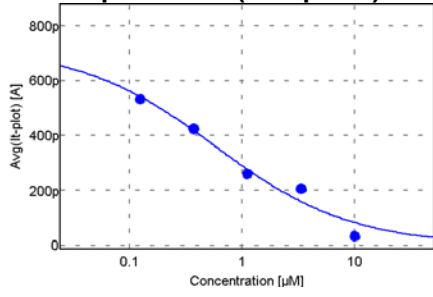
Exp 1420.4.1 (Verapamil)



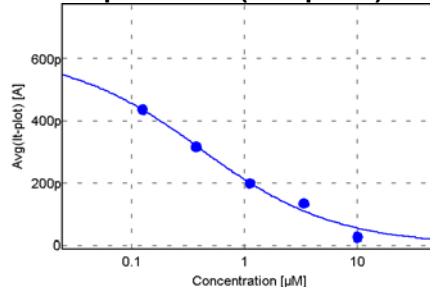
Exp 1420.4.2 (Verapamil)



**Exp 1420.4.3 (Verapamil)**

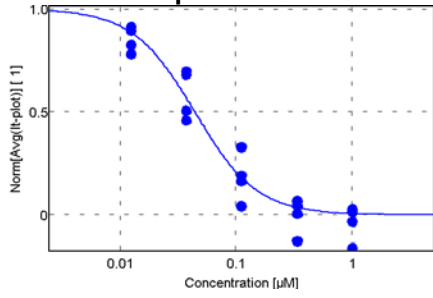


**Exp 1420.4.4 (Verapamil)**

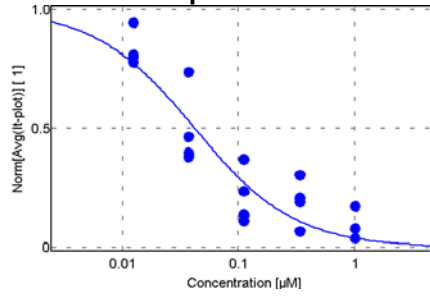


**GrpHillFit, input: Concentration, Norm[Avg(I<sub>t</sub>-plot)]**

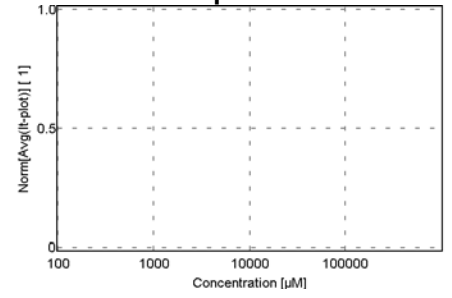
**Group Astemizole**



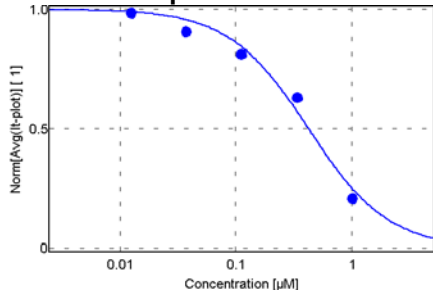
**Group Pimozide**



**Group Saline**



**Group Terfenadine**



**Group Verapamil**

