



Verification of CiPA Recommended Voltage Protocols in Patch-Clamp Assay for hERG, Cav1.2 and Late Nav1.5 Currents

Yoshimi KATAYAMA^{1),2)}, ○ Hiroshi MATSUKAWA^{1),3)}, Tomokazu KANEHISA^{1),4)}, Ayane ABE^{1),5)}, Takashi YOSHINAGA^{1),6)}, Keiichi ASAKURA^{1),7)}, Kimihito YOSHIKAWA^{1),8)}, Kazuya TSURUDOME^{1),9)}, Tadaaki TSUKAMOTO^{1),2)}

1) iSmart, Japanese Safety Pharmacology Society, 2) Pharmaceutical Research Department, Biological Research Laboratories, Nissan Chemical Corporation, 3) Higashimatsuyama Laboratories, Drug Safety Testing Center Co., Ltd., 4) Central Pharmaceutical Research Institute, Japan Tobacco Inc., 5) Drug Safety Evaluation, Research Laboratory for Development, SHIONOGI & CO., LTD., 6) Global Cardiovascular Assessment, Eisai Co., Ltd., 7) Pharmacokinetics and Safety Assessment Dept., Discovery Research Labs, NIPPON SHINYAKU CO., LTD., 8) Nonclinical Research Center, Drug Development Service Segment, LSI Medience Corporation, 9) Sophion Bioscience K.K.

Introduction

As revising an ICH regulatory guideline, S7B, for the Non-clinical Evaluation of the Potential for Delayed Ventricular Repolarization (QT Interval Prolongation by Human Pharmaceuticals) was suggested, Comprehensive *In vitro* Proarrhythmia Assay (CiPA) activities are being implemented. The suggested S7B revision includes an assay of cardiac ion channels other than hERG, and CiPA-recommended voltage protocols for patch clamping was recently announced. We, Group 5 of the JSPS iSmart (investigation of *in silico* / *in vitro* model for arrhythmogenic risk prediction) play a role of wet-data collection. This time we have investigated the applicability of those protocols.

Materials & Methods

Manual patch clamp and an automated patch-clamp system, QPatch, were selected as the platforms. The manual system was used for hERG dynamic protocol (physiological temperature) and Cav1.2 (room temperature); QPatch for hERG IC₅₀ only protocol and late Nav1.5 (room temperature). Test compounds were selected from each of the low-, intermediate-, and high-risk categories of the “training drugs” in the CiPA *in-silico* model. Flecainide was also selected.

● CiPA compounds used:

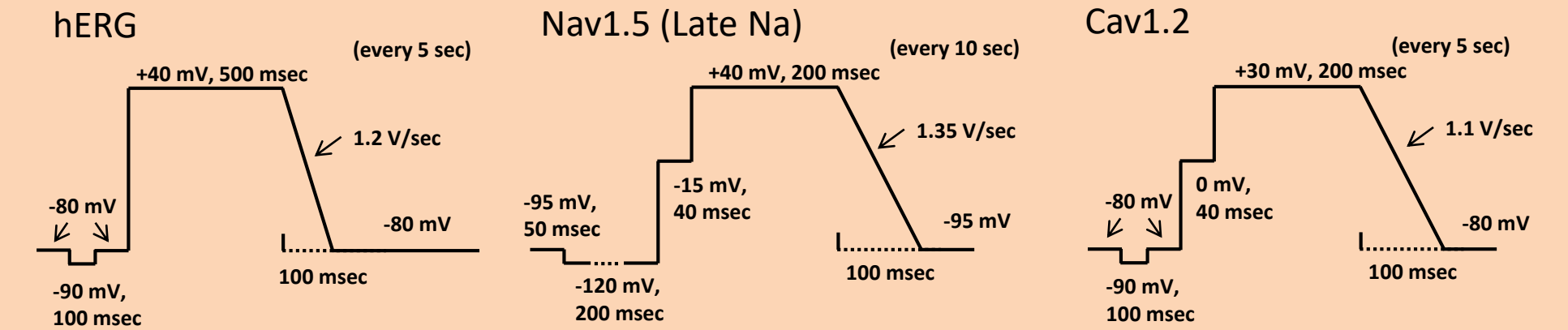
TdP risk	High	Intermediate	Low
CiPA drug	Bepidil Dofetilide	Astemizole Chlorpromazine Cisapride	Ranolazine Verapamil

 plus Flecainide

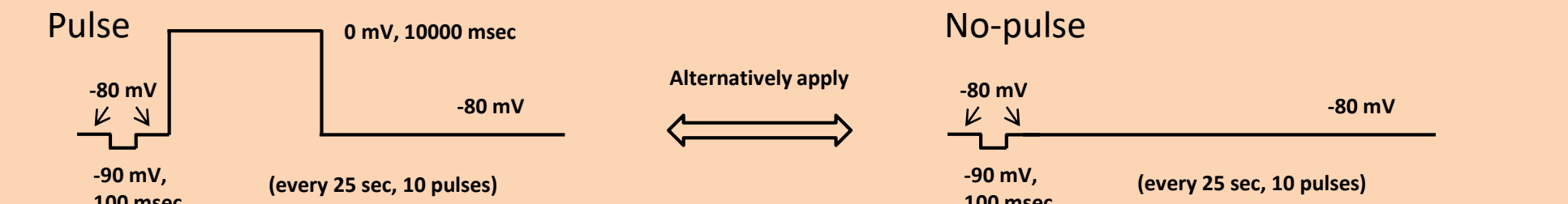
● Cell lines: hERG-HEK293/CHO-K1 cells, Nav1.5-HEK293/CHO-K1/CHL cells, Cav1.2-CHO.

● Patch-clamp conditions: identical to what the CiPA initiative announced unless otherwise stated.

➢ IC₅₀ protocols



➢ hERG dynamic model protocols



Results & Discussion

Results of hERG Assay on QPatch with IC₅₀ Only Protocol

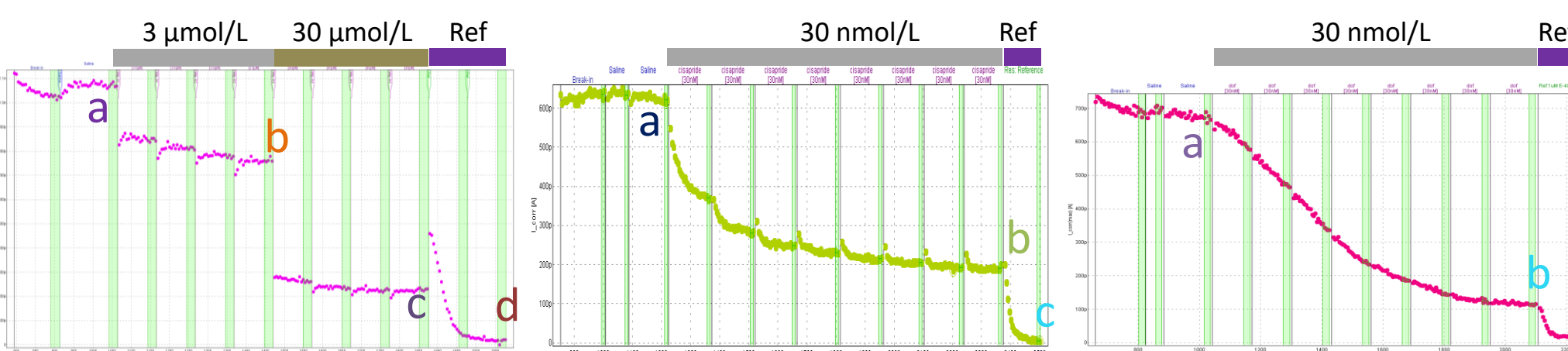
➢ I-t Plot

The compound was injected 8 times (single application) or 4 times (accumulative application) at every 20 pulses.

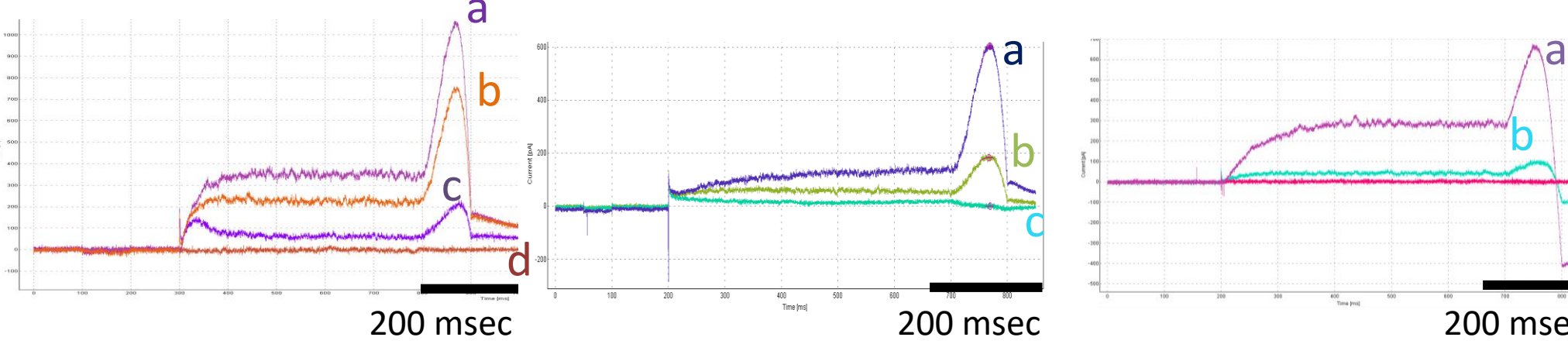
● Ranolazine

● Cisapride

● Dofetilide



➢ Current Traces



➢ Comparison with CiPA Data Set

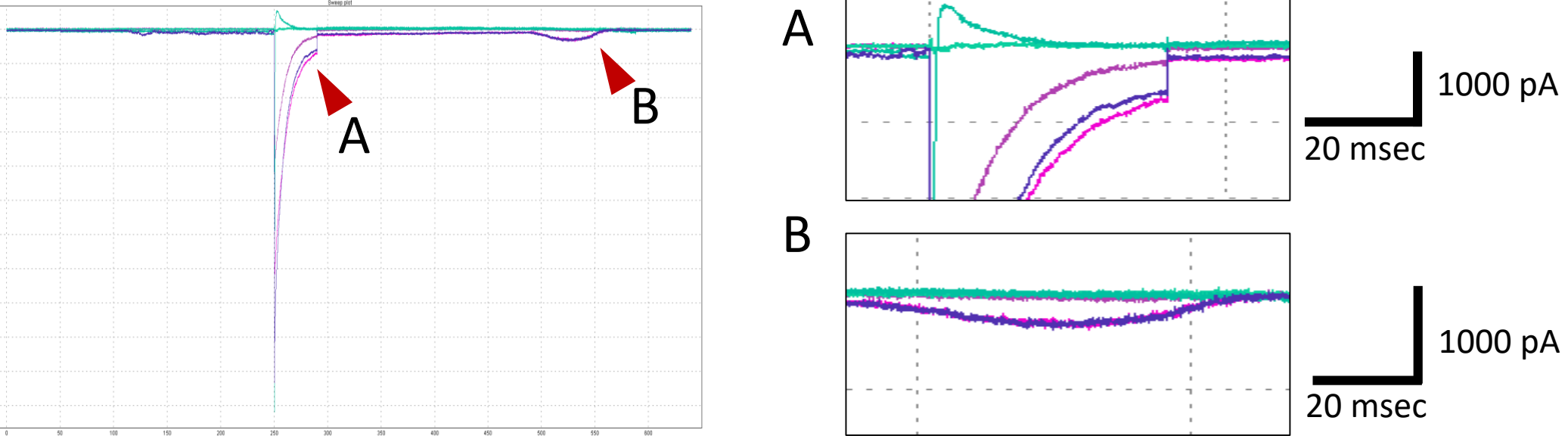
Drug	Site A			Site B		
	IC ₅₀ (μmol/L)	Hill	n	IC ₅₀ (μmol/L)	Hill	n
Bepidil	0.26	1.9	6-7	0.28	1.3	6
Dofetilide	0.015	1.5	6-8	0.011	1.8	6
Astemizole	0.026	1.4	6-9	0.026	1.3	6
Chlorpromazine	1.7	2.0	6-7	0.85	1.6	6
Cisapride	0.078	1.6	6-10	0.061	1.3	6
Ranolazine	9.0	0.92	6	9.0	0.91	6
Verapamil	0.90	2.1	7-10	0.63	1.2	6
Flecainide	1.8	1.0	7-8	1.2	0.98	6

Drug	Site C			Site D		
	IC ₅₀ (μmol/L)	Hill	n	IC ₅₀ (μmol/L)	Hill	n
Bepidil	0.031	1.3	6	0.10	1.6	7-8
Dofetilide	0.0095	1.2	6	0.014	1.8	6-8
Astemizole	0.0042	1.2	6	0.0066	1.00	6-9
Chlorpromazine	0.17	1.2	6	0.50	1.3	6-8
Cisapride	0.0080	0.93	6	0.019	0.93	6-11
Ranolazine	3.5	0.89	6	5.2	0.89	7-9
Verapamil	0.13	1.1	6	0.35	0.90	6-8
Flecainide	0.44	0.82	6	0.59	0.90	6-8

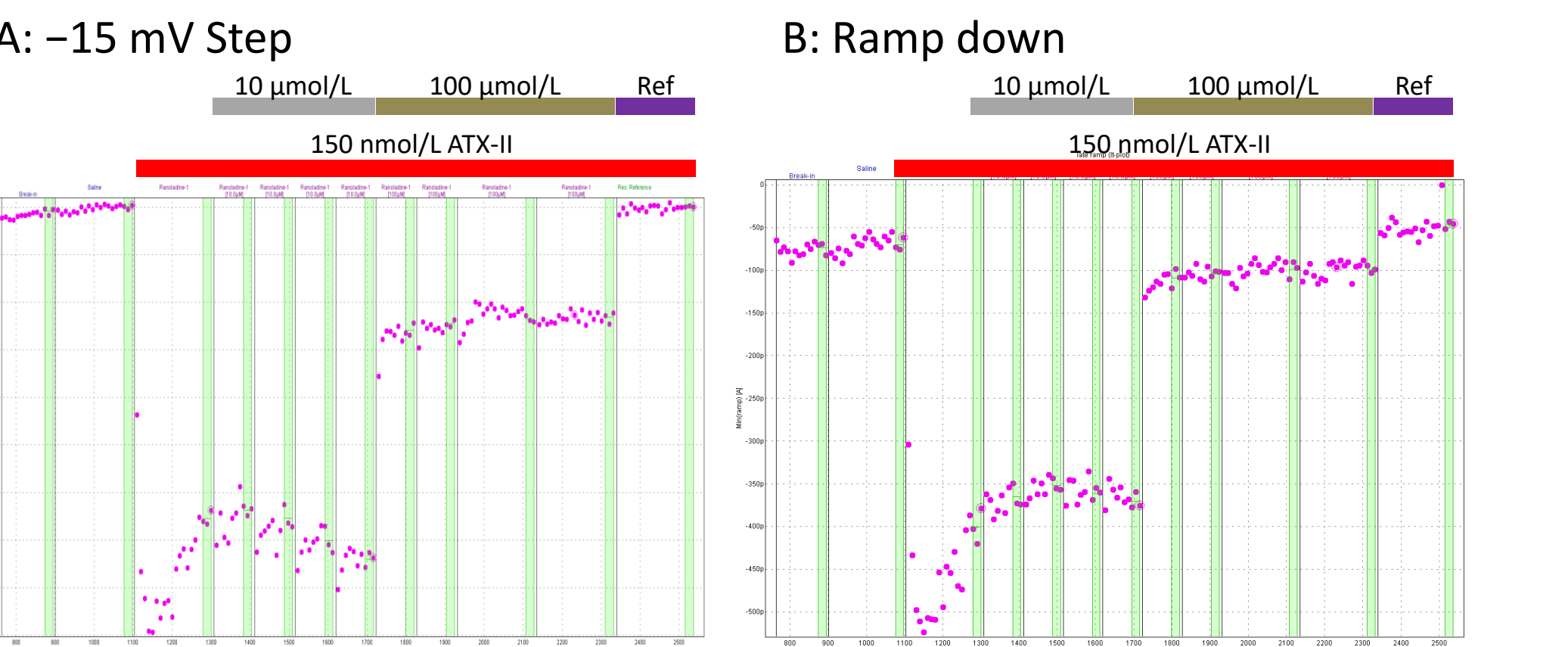
Drug	Li et al. 2018			Crumb et al. 2016		
	IC ₅₀ (μmol/L)	Hill	n	IC ₅₀ (μmol/L)	Hill	n
Bepidil	0.26	1.4	0.15	0.93		
Dofetilide	0.013	1.4	0.0015	0.63		
Astemizole	0.023	5.4	0.010	0.54		
Chlorpromazine	0.85	1.7	1.12	0.90		
Cisapride	0.071	1.9	0.012	1.3		
Ranolazine	3.4	1.1	6.5	0.84		
Verapamil	0.17	1.3	0.50	1.1		
Flecainide	—	—	—	—		

Results of Late Nav1.5 Assay on QPatch

➢ Current Traces



➢ Effect of Ranolazine on Late Nav1.5 Current

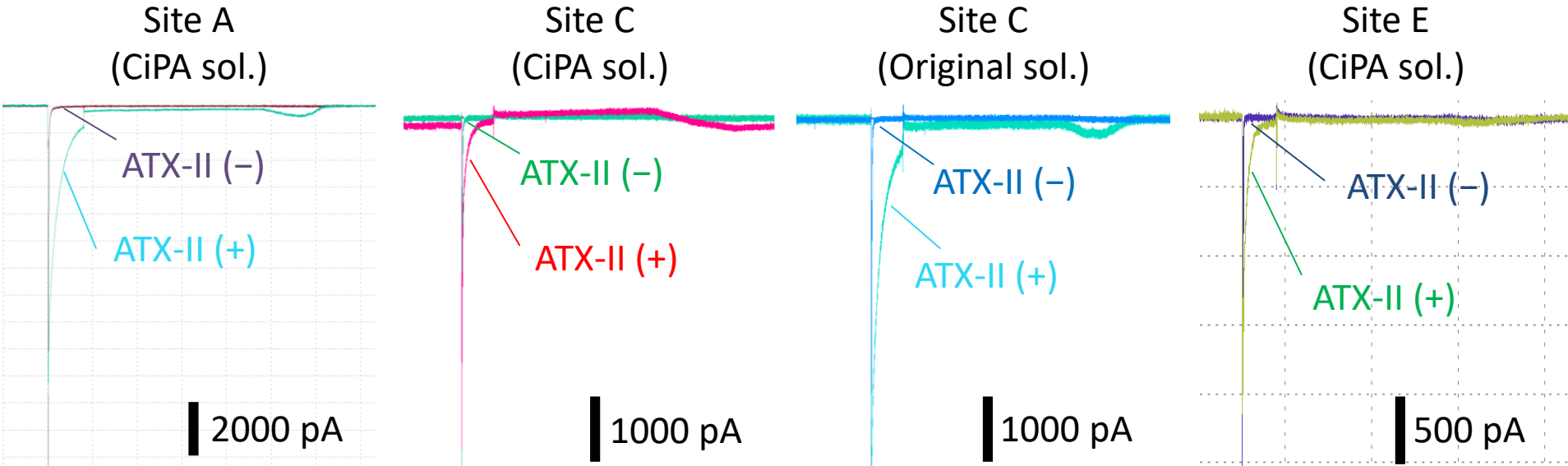


➢ Comparison with CiPA Data Set

Drug	Site A				n
	IC ₅₀ (μmol/L)	Hill	IC ₅₀ (μmol/L)	Hill	
Bepidil	1.9	1.5	1.2	0.5	3-5
Astemizole	1.7	1.8	1.1	1.3	3-6
Chlorpromazine	1.9	2.6	1.3	1.9	3-4
Ranolazine	52	1.7	29	1.7	7-8
Verapamil	14	1.4	6.3	1.3	5
Flecainide	1.5	1.4	1.3	1.0	3

Drug	Li et al. 2018		Crumb et al. 2016	
	IC ₅₀ (μmol/L)	Hill	IC ₅₀ (μmol/L)	Hill
Bepidil	0.34	1.9	1.8	1.4
Astemizole	0.60	3.1	10	2.3
Chlorpromazine	0.67	1.8	4.6	0.94
Ranolazine	6.0	0.99	7.9	0.95
Verapamil	0.98	1.2	24	2.0
Flecainide	—	—	—	—

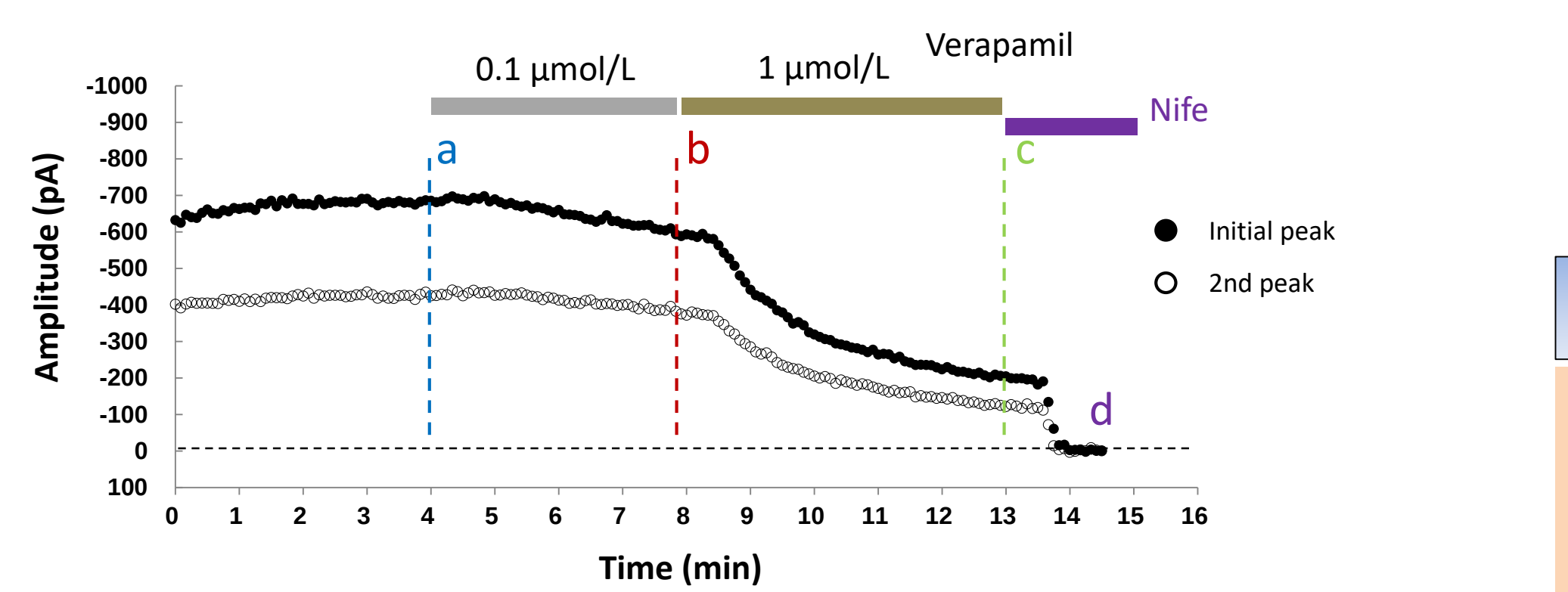
➢ Inter-facility Difference in Enhancement by ATX-II



The results show that enhancement of late Nav1.5 current by ATX-II differs between facilities and that a distinct combination of intra- or extra-cellular solutions (see below) sufficiently induced late Na current in an identical platform (site C) where late Na current could be hardly detected with CiPA conditions in the presence of ATX-II.
Note: Solutions (in mmol/L): EC, NaCl 145, KCl 4, CaCl₂ 2, MgCl₂ 1, glucose 10, HEPES 10, pH 7.4 with NaOH; IC, NaCl 10, CsF 135, EGTA 1, HEPES 10, pH 7.3 with CsOH.

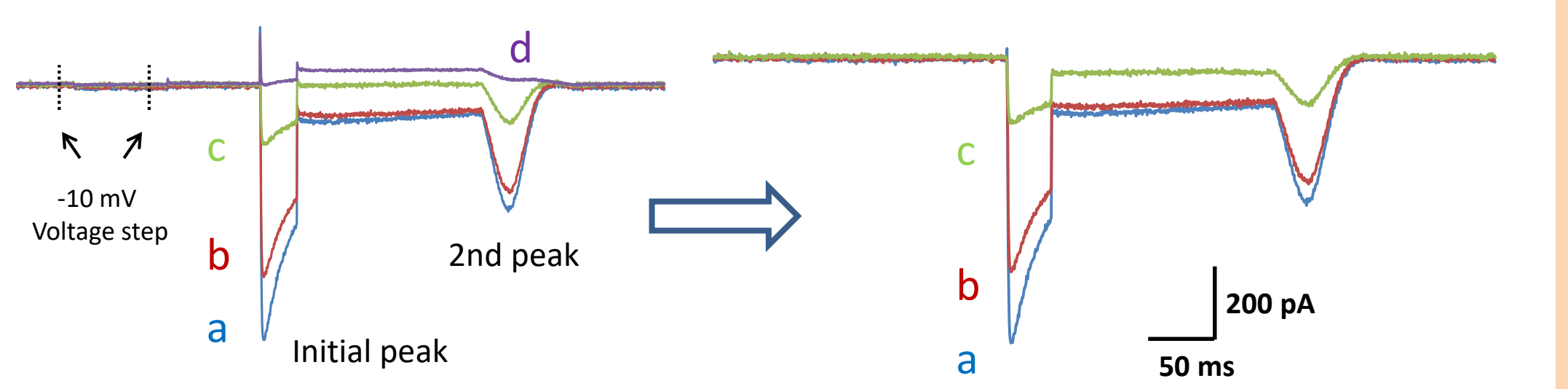
Results of Cav1.2 Assay in Manual Patch-clamp

➢ I-t Plot: Accumulative Application of Verapamil



➢ Analysis

● Cav1.2 current waveforms



● Subtraction with nifedipine treatment

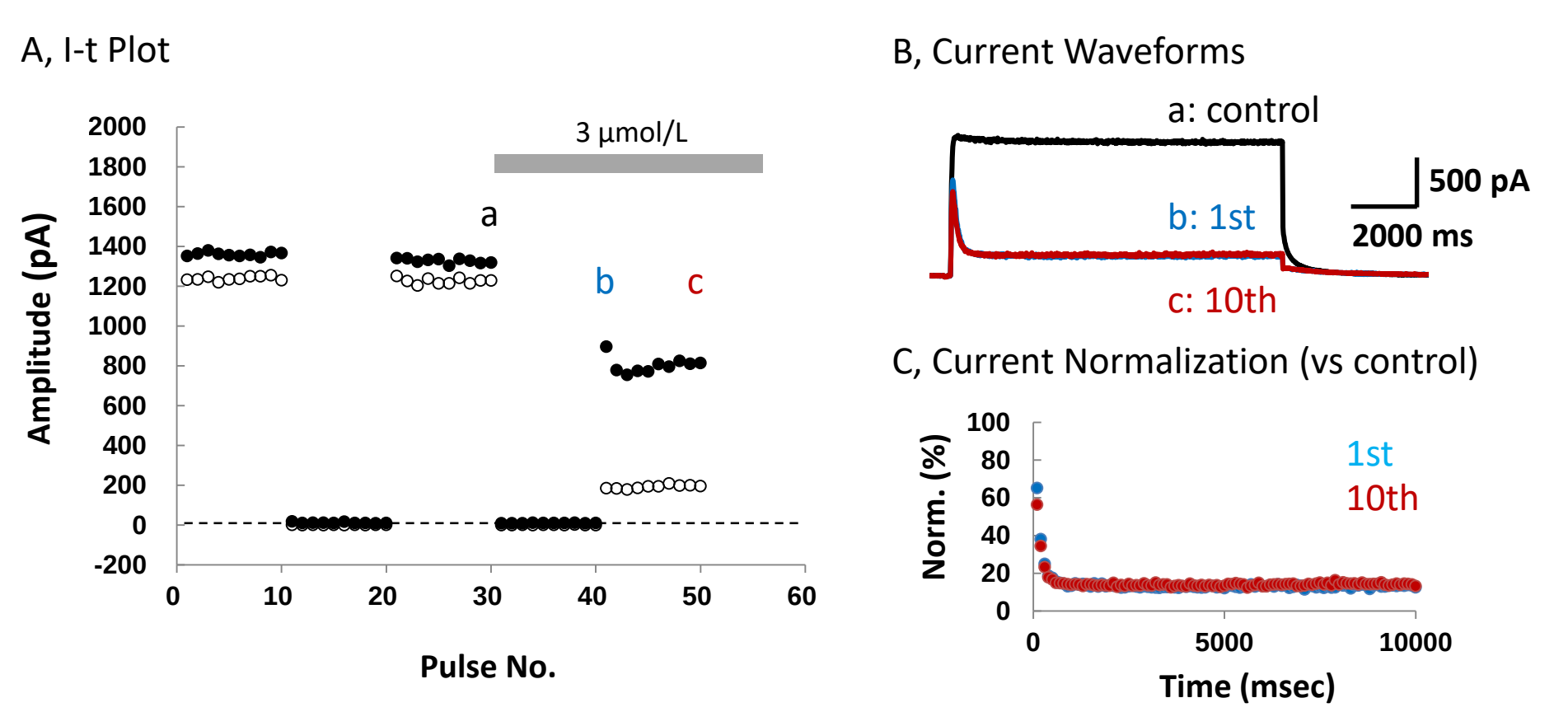
➢ Comparison with CiPA Data Set

Drug	Test conc. (μmol/L)	n	Initial peak		2nd peak		Manual ^{1,3}		HTS	
			IC ₅₀	Hill	IC ₅₀	Hill	IC ₅₀	Hill	IC ₅₀	Hill
Bepidil	0.1, 0.3, 1, 3	4	0.392	1.08	0.365	1.22	2.82	0.65	638	4.6
Dofetilide ^a	100	4	>> 100 ¹	—	>> 100 ²	—	44.5	3.6	2.30E+03	5.4
Astemizole ^a	0.03, 0.1, 0.3, 1	4	0.192	1.20	0.217	1.18	0.553	1.2	1.08	5.9
Chlorpromazine ^a	0.1, 0.3, 1, 3	4	0.741	1.59	0.728	1.98	8.32	0.85	6.35	2
Cisapride	0.3, 1, 3, 10	4	2.65	1.08	1.49	1.24	1.03E+03	4.8	4.05E+03	5.6
Ranolazine	30, 100, 300, 1000	4	514	0.99	324	0.92	900	3.9	6.54E+03	3.8
Verapamil	0.1, 0.3, 1, 3	4	0.387	0.98	0.509	1.02	0.204	1.1	11.2	0.8
Flecainide	3, 6, 30, 60	4	23.8	0.88	21.5	1.00	Crumb et al. 2016	Li et al. 2018		

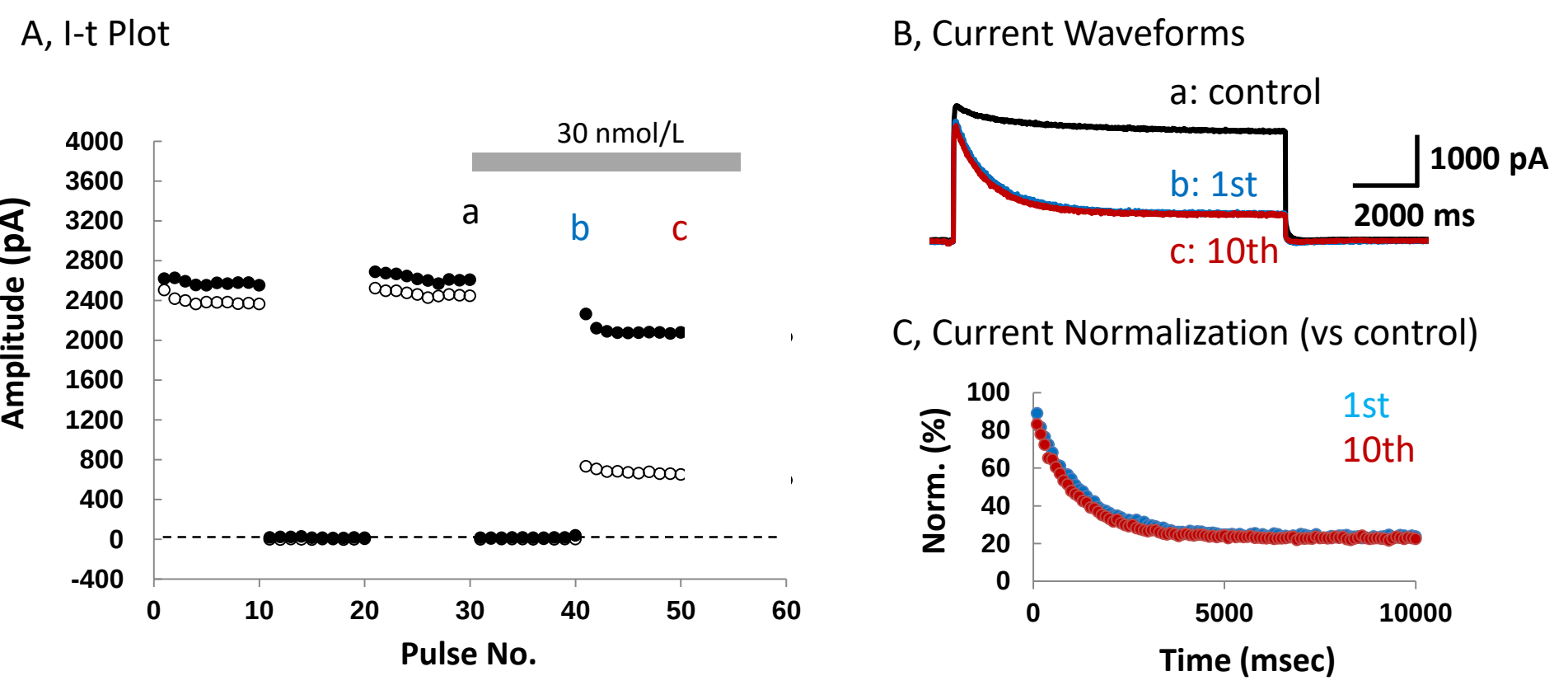
*1: 8.6% at 100 μmol/L. *2: 5.7% at 100 μmol/L. *3: Action potential protocol, Ba²⁺ as charge carrier, at physiological temperature.
#: Each cell was exposed to single drug concentration.

Results of hERG Dynamic Protocol

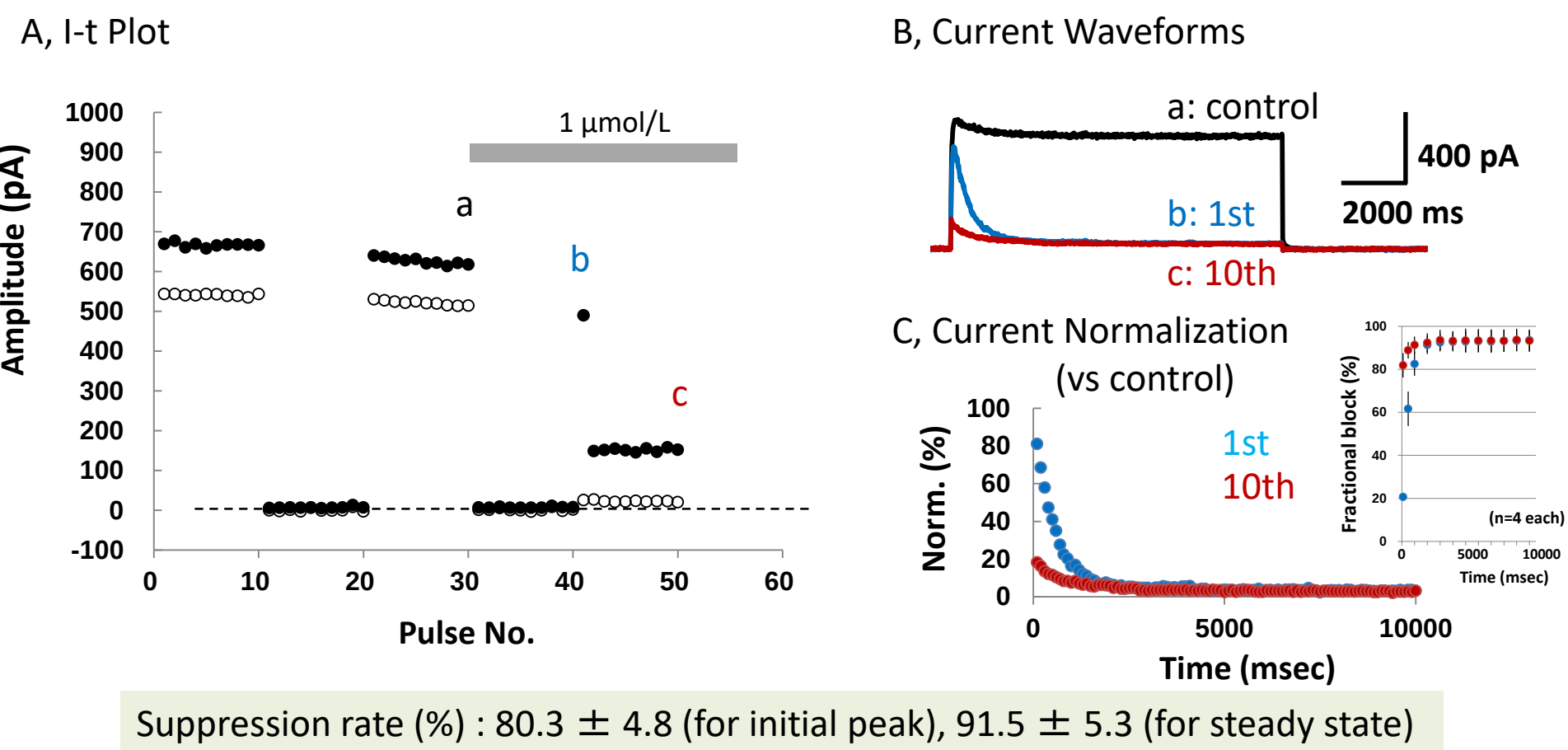
➢ Effect of Verapamil on hERG Current



➢ Effect of Cisapride on hERG Current

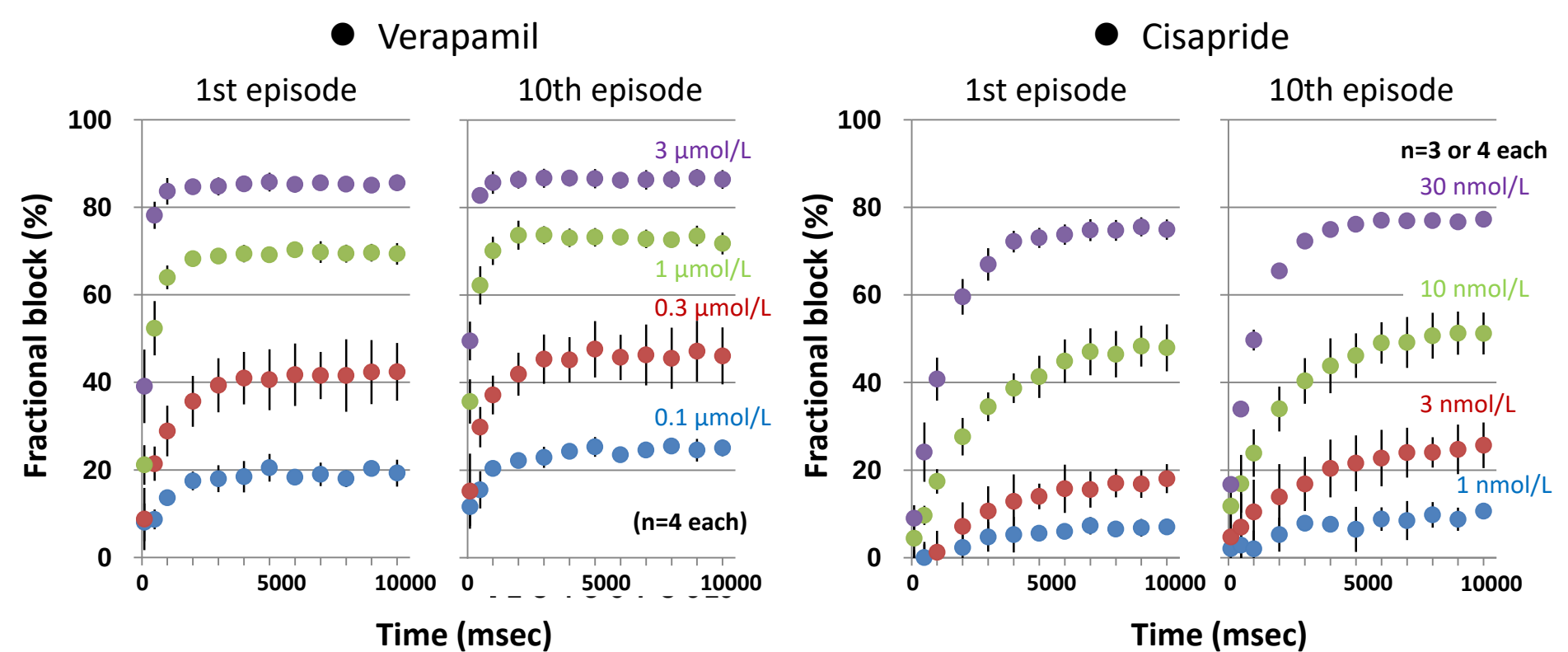


➢ Effect of E-4031 on hERG Current



Suppression rate (%) : 80.3 ± 4.8 (for initial peak), 91.5 ± 5.3 (for steady state)

➢ Concentration-dependent Progressive Block



➢ Comparison with CiPA Data Set

Drug	Test concentration (μmol/L)	n	Initial peak IC ₅₀ (μmol/L)	Hill	Steady state IC ₅₀ (μmol/L)	Hill	IC ₅₀ (μmol/L)	Hill
Cisapride	1, 3, 10, 30*	3 or 4	NA	NA	10.8 ²	1.11	10.1 ¹	0.7
Ranolazine	1, 3, 10, 30	3 or 4	13.5 ¹	0.67	8.50	0.89	8.27 ¹	0.9
Verapamil	0.1, 0.3, 1, 3	4	2.48	0.60	0.369 ¹	0.85	0.288 ¹	1
Flecainide	0.3, 1, 3, 10	4	2.47	0.65	1.16	0.86	Li et al. 2017, 2018	

*: Cisapride concentrations are shown in nmol/L. NA: Not applicable.

Conclusions

- Some of our data which disagree with those presented by CiPA are discussed. Some problem and suggested modifications are pointed. We plan to deliver these suggestions to CiPA.
- In hERG IC₅₀ only protocol, the similar results were obtained among different facilities after optimizing the duration and number of application on QPatch system.
- In late Nav1.5 current protocol, it was turned out that enhancement of late Na current by ATX-II was not consistent among different facilities when CiPA solutions were used and was dependent on the combination of intra- and extra-cellular solutions, suggesting the possibility that the protocol may still have some room for improvement.
- In Cav1.2 current protocol, the IC₅₀ values of dofetilide and cisapride deviated far from the results by CiPA.
- In a limited number of compounds from intermediate and low risk categories, the IC₅₀ values were almost similar to the results by CiPA.

References

- Recommended voltage protocols to study drug-cardiac ion channel interactions using recombinant cell lines
- Journal of Pharmacological and Toxicological Methods, Crunb et al., 2016
- Circ Arrhythm Electrophysiol, Li et al., 2017
- CLINICAL PHARMACOLOGY & THERAPEUTICS, Li et al., 2018

Acknowledgement

We would like to express our gratitude to Sophion Bioscience K.K. for the courtesy of QPlates and the solid technical support with respect to data acquisition and analyses on QPatch.