

Introduction

The voltage dependent sodium channel is responsible for the upstroke and directed propagation of action potentials in nerve and muscle cells and is therefore a central ion channel in excitable tissues. The implication of voltage gated sodium channels in pain mediation and diseases such as epilepsy and cardiac arrhythmia has made them very important targets for drug discovery. Here we show eight subtypes of the voltage gated sodium channel tested in parallel on Qube, the Sophion 384format automated patch clamp system designed for high throughput and high fidelity electrophysiological recordings. On Qube it is possible to run up to 16 different clones or cell lines simultaneously which ensures identical conditions across the experiments.

Three types of experiments were designed to explore 1) TTX sensitivity, 2) IV-relationship for activation and inactivation and 3) pulse train suitable for screening for use dependent sodium channels blockers. For the screening example pharmacology was represented by tetracaine. For every run Na_v channel subtypes Na_v1.1, 1.2, 1.3, 1.4, 1.5, 1.6, 1.7 and 1.8 were tested side by side.



Na _v 1.1 -	С ^А в	<u> </u>	ار ا		1	ייר קייי	Т 7	1	**	1	9 9	"		T T	T	"" "	7	~	1	7	***	┓	***		***
Na _v 1.2 -		T	7-	T 7~	7	77	1	***	1	1	7		1 -	<u>ד</u>	T	***	***		77	1	7		**	7	
Na _v 1.3 -	E F	1	<u>-</u>	T	- -	4	1		- -	**				1		<u>۳</u>	1			**	**	<u>~</u>		<u>1</u>	7
Na _v 1.4 -	С ч	É.				1 1 1		<u> </u>	- - -			1 1 1	77	2	77	- 					1		T		T
Na _v 1.5 -				7		<u> </u>						****											<u> </u>	7 7	
• Na√1.6 ⊀	к												7 7 7											<u>r</u>	1
• Na√1.7 ⊰				1												-									
Na. 1.8		7	, , ,		7	Y	Y	r	r	r		1	7	7	1 	r	r		۲ ۲	7		r	7	r	1
	L P		r	~	Y	Y	Y	r	1	۳ ۲	r	Υ-	Y	r	r	1	Υ	r	Y	r	m	r	Υ	Y	~ ~

1 2 3 4 5 6 7 8 9 10 11 12 13 14 15 16 17 18 19 20 21 22 23 24

[Fig. 1]: Top, cell clone Cell Transfer Plate (ccCTP) for the Qube instrument. Cells expressing the different sodium channel isoforms were added to the individual wells of the ccCTP and subsequently transferred to the QChip for experiment. Lower panel shows layout and evoked currents from a typical experiment. Currents were elicited by a depolarizing pulse from a holding potential of -110 mV to 0 mV.



References

Sophion Qube

Catteral W.A. et al. Pharmacol Rev. 57:397-409, 2005 and references therein.







response to increasing doses of TTX.



Materials and methods

Cells expressing sodium channel isoforms Nav1.1 to Nav1.8 were cultured according to the SOP for the respective Solutions: Intracellular solution (IC) (in mM): CsF 135, NaCl 10, EGTA 5, HEPES 10, adjusted to pH 7.2 (CsOH). cell line. Nav1.1, Nav1.2, Nav1.3 and Nav1.4 were from SB Drug Discovery (Glasgow, UK), Nav1.5 was from B'SYS Extracellular solution (EC) (in mM): NaCl 145, KCl 4, MgCl₂ 1, CaCl₂ 2, HEPES 10, glucose 10, adjusted to pH 7.4 GmbH (Witterswil, CH), Nav1.6 and Nav1.8 from ChanTest (Cleveland, OH) and Nav1.7 from Anaxon AG (Berne, (NaOH). CH). On the day of experiment the cells were harvested using detachin or trypsin and transferred to serum-free medium (EX-CELL® ACF CHO Medium, Sigma-Aldrich, Brøndby, DK) supplemented with HEPES 25 mM, 40 µg/ml trypsin inhibitor and P/S. At the time of the experiment, the cells were washed with EC solution and placed in the Analysis was carried out using the Sophion Analyzer software and the R package "drc"x (www.R-project.org and cell clone Cell Transfer Plate. www.bioassay.dk).



Biophysical and pharmacological characterisation of multiple Na_v subtypes on Qube

Anders Lindqvist, Søren Friis, Mette T. Christensen & Mads P.G. Korsgaard

[Fig. 2]: Voltage-dependent activation and steady-state inactivation for Na_v1.X isoforms. Figure 2A, typical currents in response to a depolarizing pulse from -110 to 0 mV and figure 2D, voltage protocol. Figure 2B, conductance-voltage relationship for activation and steady-state inactivation for Na_v1.2 (blue), Na_v1.6 (orange) and Na_v1.7 (grey). Figure 2C, activation and steady-state in-activation for Na_v1.7 under control conditions (black) and in the presence of 10 μM of amitriptyline (blue), bupivacaine (orange) and tetracaine (grey) respectively.



[Fig. 4]: Dose-response relationship in response to tetracaine. Fig. 4A, whole-cell currents in the presence of increasing concentrations of tetracaine. Insert shows details of currents from Nav1.5 and Nav1.6 expressing cells. Figure 4B, voltage protocol. Currents were measured at first and tenth (last) depolarization. Figure 4C, dose-response relationships for Nav1.1 (blue), Nav1.4 (orange)

Conclusion

This study presented biophysical and pharmacological characterization of eight different Nav cell lines on Qube simultaneously. Multiple compounds and cell lines were used in parallel on Qube and the results for Tetrodotoxin sensitivity, IV-relationship (activation and inactivation), and screening of use dependent sodium channel blockers were of high and consistent quality, which proves excellent voltage control and cell handling. This Qube feature creates endless opportunities for simultaneous testing of multiple cell lines and compounds on Qube.

[Table 1]: Half maximal activation and inactivation for the respective channel isoform.

Channel isoform	V½ activation mean ± SEM (mV)	V½ inactivation mean ± SEM (mV)	V½ inactivation at 10 μM tetracaine (mV)	n
Na _v 1.1	-29 ± 0.4	-56 ± 0.7	-77 ± 1.2	10
Na _v 1.2	-26 ± 0.4	-56± 0.4	-76 ± 2.0	8
Na _v 1.3	-30± 0.3	-66± 0.4	-85 ± 0.5	6 - 8
Na _v 1.4	-35± 0.4	-67± 0.4	-84 ± 0.2	3
Na _v 1.5	-36± 1	-60± 0.9	-77 ± 1.6	12
Na _v 1.6	-36± 0.5	-63± 0.3	-85 ± 0.3	8
Na _v 1.7	-36± 0.4	-66± 0.2	-86 ± 0.2	10
Na _v 1.8	-11± 0.3	-45± 0.4	-63 ± 1.3	10

[Table 2]: Concentration dependent inhibition of sodium currents by TTX. Cells were stimulated with depolarizations from -110 to 0 mV at 0.1 Hz.

Channel isoform	IC ₅₀ ± SEM (nM)	n
Na _v 1.1	22 ± 5	11 - 14
Na _v 1.2	33 ± 4	11 - 15
Na _v 1.3	12 ± 3	7 -14
Na _v 1.4	44 ± 5	7 - 13
Na _v 1.5	7800 ±	7 - 12
Na _v 1.6	9 ± 1	11 - 15
Na _v 1.7	93 ± 9	12 - 15
Na _v 1.8	>10000	13 - 16

[Table 3]: Concentration-dependent inhibition of sodium currents by tetracaine. Cells were stimulated with 10 depolarizations from -110 mV to 0 mV at 10 Hz.

Channel isoform	First depolarization $IC_{50} \pm S.E.M (\mu M)$	Tenth depolarization IC ₅₀ ± S.E.M (µM)*	n
Na _v 1.1	>30	8.6 ± 2.8	2 - 8
Na _v 1.2	>30	9.7 ± 2.9	4 - 6
Na _v 1.3	>30	9.9 ± 0.3	2 - 8
Na _v 1.4	>30	3.4 ± 0.4	2 - 4
Na _v 1.5	>30	3.1± 0.2	5 - 9
Na _v 1.6	>30	11.4 ± 4.9	6 - 8
Na _v 1.7	>30	3.9 ± 0.6	3 - 7
Na _v 1.8	>30	4.3 ± 3.4	4 - 7