

ENDOGENOUS ION CHANNELS OF MAMMALIAN CELL LINES CHARACTERIZED WITH THE QPATCH™



SOPHION BIOSCIENCE A/S
Baltorpevej 154
DK-2750 Ballerup
DENMARK
info@sophion.dk
www.sophion.com

SOPHION BIOSCIENCE, INC.
675 US Highway One
North Brunswick, NJ 08902
USA
Phone: +1 732 745 0221
www.sophion.com

¹SØREN FRIIS

¹RASMUS B. JACOBSEN

¹MORTEN R. SUNESEN

¹DORTHE NIELSEN

¹NADIA LARSEN

²CHRIS MATHES

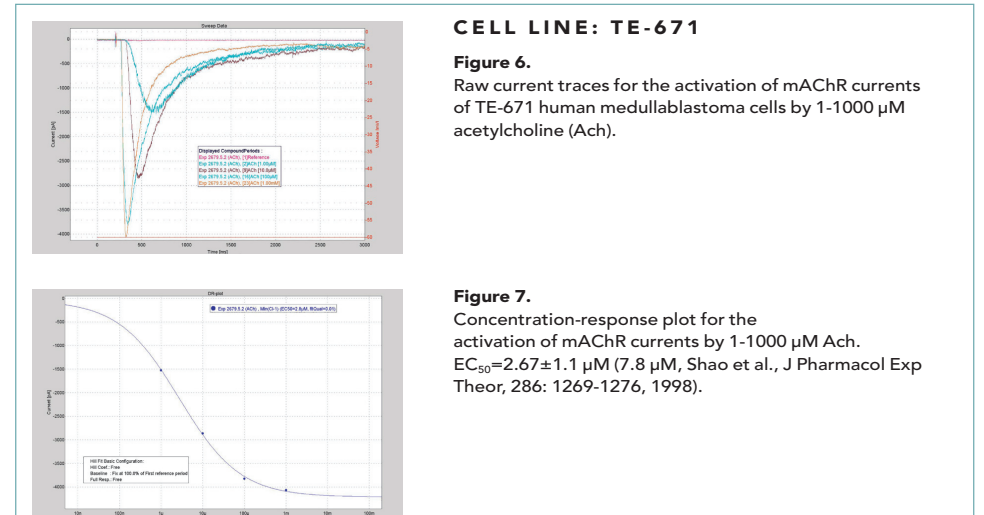
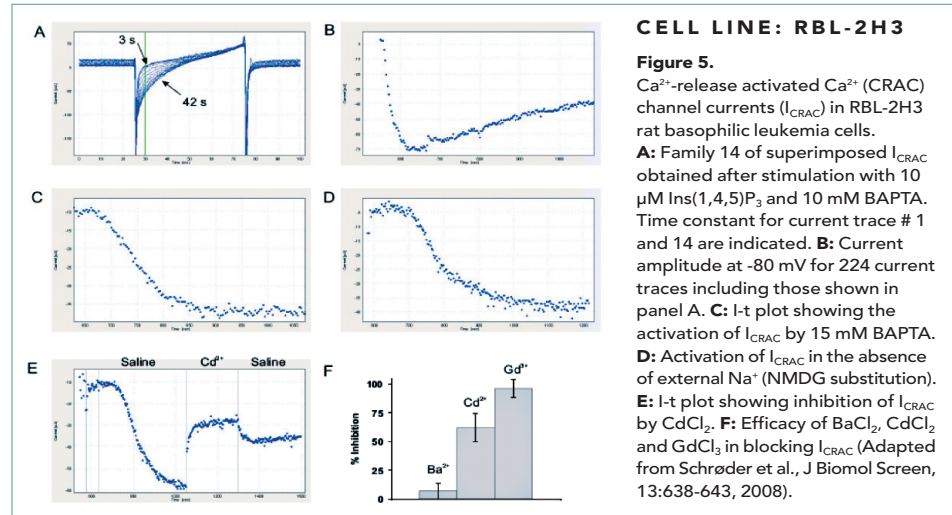
³NIELS J. WILLUMSEN

¹Sophion Bioscience A/S, 154 Baltorpevej, DK-2750 Ballerup, Denmark. | ²Sophion Bioscience, Inc., North Brunswick, NJ 08902, USA | ³Department of Biology, University of Copenhagen, DK-2100 Ø, Denmark

A large number of mammalian cell lines are commercially available to be used as expression systems for membrane or cytoplasmic proteins. A number of voltage and ligand gated ion channels of potential interest for the pharmaceutical industry are endogenously expressed in several CNS and non-CNS cell lines including TTX-sensitive Na⁺ channels, Ca²⁺-release activated Ca²⁺ (CRAC) channels, inward rectifier K⁺ channels, acid-sensing ion channels (ASIC) and muscarinic alpha-adrenergic receptors mAChR). We have explored the applicability of five commonly employed cell lines from American Type Culture Collection (ATCC) for use with Sophions QPatch™ automated patch clamp systems (QPatch 16 and QPatch HT) and characterized the ion channel types that they endogenously express. Specifically we have explored:

- Suitability for automated patch clamp studies ('patchability')
- Background ionic currents that may interfere with currents of experimentally expressed ion channels
- Possible use for characterizing ion channels of interest without the need to experimentally introduce their genes (expression)

The tests have led to the development of a number of simple standard operation procedures (SOPs) for employment of the cell lines in QPatch characterizations of ion channels.



BASIC PROPERTIES OF FIVE CELL LINES: PATCHABILITY AND EXPRESSION OF ENDOGENOUS ION CHANNELS

Cell line	Origin	ATCC #	Basic patchability			Na ⁺ current		K ⁺ current		Ca ²⁺ current		LGIC current	
			R _{seal} (GΩ)	R _{wholecell} (GΩ)	Wholecells (%)	Expression (%)	I _{peak} (pA)	Expression (%)	I _{peak} (pA)	Expression (%)	I _{peak} (pA)	Expression (%)	I _{peak} (pA)
SH-SY5Y	Human neuroblastoma	CRL-2266™	3.42±1.01	2.08±1.01	38	28 (TTX insensitive)	-409±101	16 (K _v 3.x)	1040±358	3	-180	-	-
PC-12	Rat adrenal gland	CRL-1721™	1.22±0.35	6.47±2.30	46	-	-	53	263±164	-	-	-	-
IMR-32	Human neuroblastoma	CCL-127™	6.36±2.26	4.31±1.75	33	74	-377±68	21	113±13	-	-	-	-
RBL-2H3	Rat basophilic leukemia	CRL-2256™	1.60±0.21	1.98±0.46	77	-	-	100 K _{ir}	-1087±79 ¹	95 CRAC	-32±3	-	-
TE671	Human medullablastoma	CRL-8805™	0.87±0.16	0.87±0.13	85	100	1525±330	-	-	-	-	100 mAChR ²	1190±149

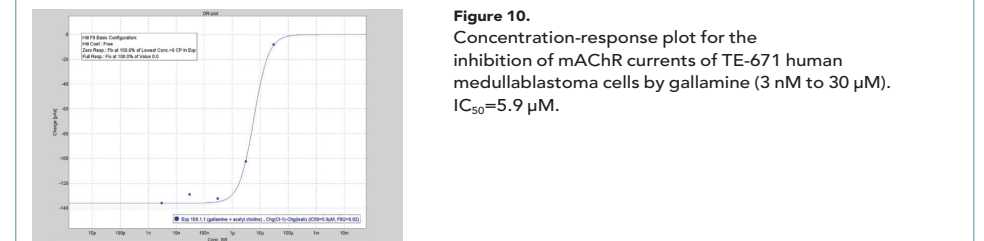
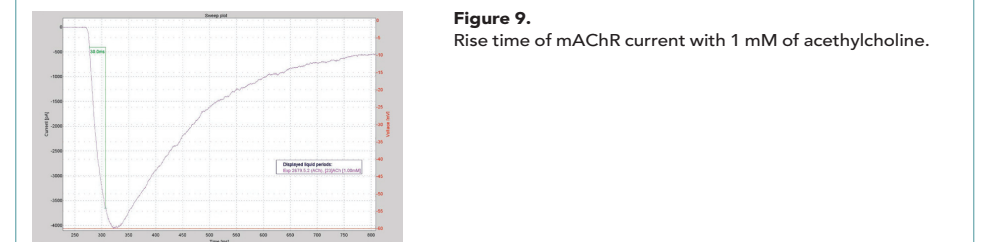
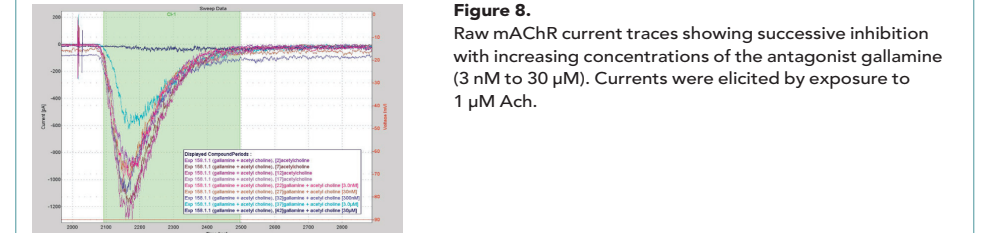
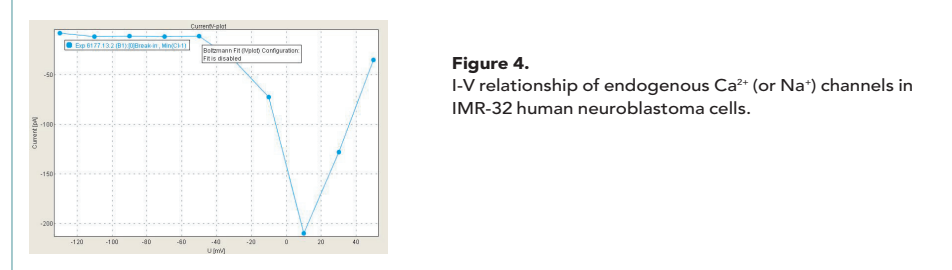
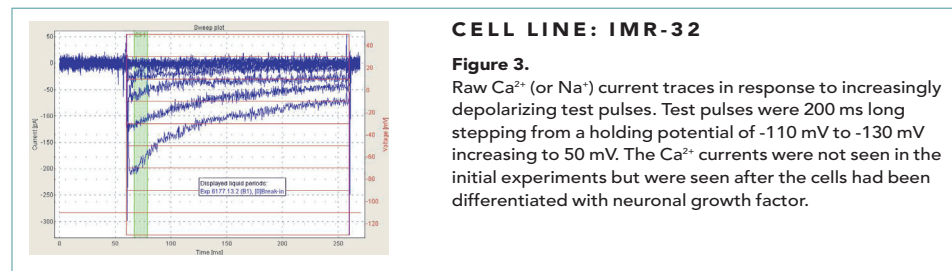
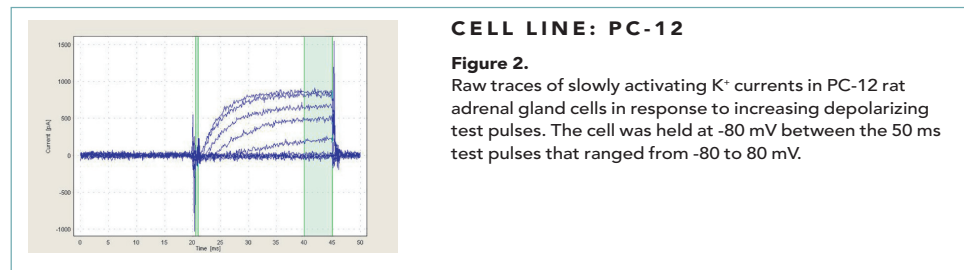
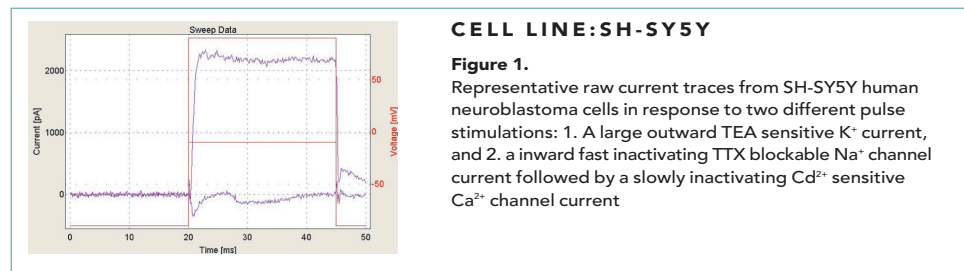
Basic Properties

The table collects the following types of data for the five cell lines that were characterized:

1. patchability with QPatch™ based on gigaseal and whole-cell resistances and whole-cell success rate
2. expression level of the endogenous ion channel species as percentage of cell recordings that exhibited the specific ionic current
3. whole-cell current amplitudes of the identified ion channels. The nature of the ion channels were based on use of blockers, activators and in substitutions. The functional properties are presented in the figures below

The patchability was tested with physiological Ringer's solutions and without specific optimization of test conditions. For all cell lines: n>40. Where the identity of ion channels is positively known the name is given without parenthesis. Ion channels listed in parenthesis indicate most likely identity.

- 1: In response to 140 mM K⁺ at -40 mV
- 2: In response to 370 μM ACh



SUMMARY:

Five commercially available cell lines were tested for use with QPatch™ automated patch clamp systems and for their expression of endogenous ion channels. All cell lines proved to be well suited for exploration with the QPatch technology. It was found that each cell line exhibited a specific and characteristic expression of ion channels including sodium, potassium and calcium channels as well as ligand-gated cation channels. It is concluded that the selection of a proper cell line in several cases may eliminate the need for expression of specific ion channel genes in drug testing assays. Standard SOPs for assays based on QPatch™ recordings of ion channel currents of the five tested, commercially available cell lines are available from Sophion Bioscience.