

Introduction

Automated patch clamp platforms have become an indispensable tool for increasing throughput for hit detection and compound validation in drug screening. As demonstrated by Turner et al. 2016, QPatch is a useful tool in characterizing CFTR functionality.

Here we demonstrate how QPatch HTX can be used for evaluation of pharmacological effects. QPatch HTX is a 48 channel format automated electrophysiology system with state-of-the-art microfluidic channels system in the QPlate.

The Cystic Fibrosis Transmembrane Conductance Regulator (CFTR) is a chloride ion channel that conducts Cl⁻ ions across epithelial cell membranes. The CFTR channel regulates fluid movement and is regulated by phosphorylation. The CFTR channel consists of two transmembrane domain regions and has three intracellular domains that regulate channel activity, two nucleotide binding domains and the R-domain (Regulatory domain). The gating of the CFTR channel requires both phosphorylation of PKA (cAMP-dependent protein kinase) at the unique Regulatory domain and ATP binding and hydrolysis at the Nucleotide binding site (Carson et al. 1995).

Forskolin is an activator of the CFTR channel. Forskolin works via activation of adenylate cyclase that causes ATP to generate cAMP, which in turn works on the cAMP-dependent protein kinase (PKA). PKA phosphorylates the CFTR channel and thereby opens the channel (Haws et al. 2002). Genestein is then added to keep the channel open.

Berger et al 1998, has described how fluoride (F⁻) can stimulate the CFTR channel activity. Fluoride interferes through the adenylate cyclase signaling pathway, but by a different entry point than Forskolin. When PKA phosphorylates the CFTR channel with onset by F⁻ the channel stays open (Berger et al. 1998).

CFTR channels are constitutively open once they have been exposed to reagents activating the adenylate cyclase, which in these experiments were F⁻. Therefore is was not possible to compensate for C_{fast}, C_{slow}, R_{serie} and determinate R_{membrane} Instead a reference sweep recorded at the end of each experiment was subtracted all other sweeps in the same experiment.

Materials and methods

The cell line CHO-hCFTR was kindly supplied by ChanTest – part of Charles River Laboratories International, Inc.

The cells were cultured according to the Sophion standard operating procedure for CHO cells and harvested using Detachin and kept in the QStirrer on the QPatch prior to the automatic preparation.

Solutions

Intracellular saline solution: 2.7 mM CaCl₂, 0.88 mM MgCl₂, 10 mM HEPES, 60 mM KF, 70 mM KCl, 15.6 mM KOH/10 mM EGTA, 2 mM Na₂-ATP, pH = 7.2, 295 mOsm.

Extracellular saline solution:

1 mM CaCl₂, 1 mM MgCl₂, 5 mM HEPES, 5 mM MES, 3mM KCl, 140 mM NaCl, 20 mM TEA-Cl, pH = 7.3, 290 mOsm.

Seal- and whole cell formation

After QPatch had transferred the cell suspension to the QPlate, a pressure -70 mBar was applied to obtain positioning and sealing of the cells. A whole cell protocol with pressure pulses at -150 mBar was used to obtain whole cell formation. Thereafter the membrane was clamped at -90 mV until and between the voltage protocols were executed.

Conclusion

These experiments demonstrate that reliable recordings with the chloride conducting ion channel CFTR can be performed on QPatch using fluoride as activator of adenylate cyclase.

We have shown the ability to find the IC₅₀ for CFTR_{inh}-172 in close correspondence to the literature values.

We can conclude that the overall success in obtaining stable seals and completed experiments are at a level that clearly identifies the CHO–hCFTR assay as feasible and with high quality results on QPatch.



CFTR pharmacology determined with QPatch

Naja M. Sørensen, Mads P. G. Korsgaard, Rasmus B. Jacobsen





[Fig. 2]: Current traces during baseline conditions. The Blue arrow indicates $E_{rev} = -20 \text{ mV}$

Five point dose response using CFTR-inh 172 on CFTR



[Fig. 3]: Raw data traces in the absence as well as presence of increasing concentrations of CFTR_{inh}-172



[Fig. 5]: Hill fit for dose response with CFTR_{inh}-172, n=18. IC₅₀ = 1.55 μ M ± 0.19. Error bars are SEM

References

Carson MR, Welsh MJ. Structural and functional similarities between the nucleotide-binding domains of CFTR and GTP-binding proteins. Biophys J. (1995) ;69(6):2443-8. Haws CM, Nepomuceno IB, Krouse ME, Wakelee H, Law T, Xia Y, Nguyen H, Wine JJ. Delta F508-CFTR channels: kinetics, activation by forskolin, and potentiation by xanthines. The Journal of Biological Chemistry, October 4, (2002), 277, 37235-37241.

Berger HA, Travis SM, Welsh MJ. Fluoride stimulates cystic fibrosis transmembrane conductance regulator Cl⁻ channel activity. Am J Physiol. (1998);274(3 Pt 1):L305-12. Tonghui Ma, Jay R. Thiagarajah, Hong Yang, Nitin D. Sonawane, Chiara Folli, Luis J.V. Galietta, and A.S. Verkman. Thiazolidinone CFTR inhibitor identified by high-throughput screening blocks cholera toxin-induced intestinal fluid secretion. J Clin Invest. (2002); 110(11): 1651–1658.

Taddei A, Folli C, Zegarra-Moran O, Fanen P, Verkman AS, Galietta LJV. Altered channel gating mechanism for CFTR inhibeted by a high-affinity thiazolidinone blocker. FEBS letters 558 (2004) 52-56. Mark J. Turner, Elizabeth Matthes, Arnaud Billet, Amy J. Ferguson, David Y. Thomas, Scott H. Randell, Lawrence E. Ostrowski, Kathy Abbott-Banner and John W. Hanrahan. The dual phosphodiesterase 3 and 4 inhibitor RPL554 stimulates CFTR and ciliary beating in primary cultures of bronchial epithelia. Am J Physiol Lung Cell Mol Physiol 310: L59-L70, 2016.