

# Application Report

## RD(TE671)nAChR $\alpha 1$ on QPatch

Endogenous nicotinic channels

### Summary

The human cell line RD(TE671) endogenously expresses the nicotinic acetylcholine  $\alpha 1$  receptor (nAChR $\alpha 1$ ). The QPatch can reliably measure ligand-gated RD(TE671) currents and was used to confirm the specific activation of nAChR $\alpha 1$  due to acetylcholine addition on the QPatch.

- Stable seals rates above 80% were repeatedly obtained
- QPatch performed reliable dose-response experiments on the fast ligand-gated channel nAChR $\alpha 1$
- It was confirmed that the observed currents indeed are due to the activation of acetylcholine receptors

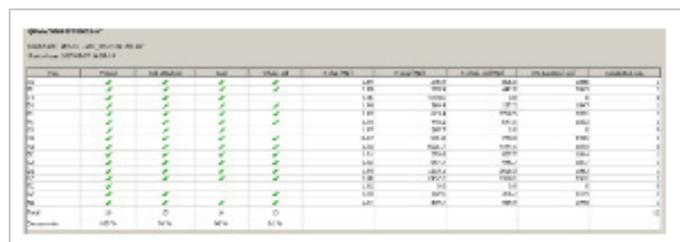


Fig. 1: Typical overview of success rates with the RD(TE671) cell line.

### Results and discussion

The RD(TE671) cells performed well on QPatch with a seal rate constantly above 80% (Figure 1). Cells were homogenous in size around 15 pF and all exhibited significant nAChR $\alpha 1$  current upon stimulation with acetylcholine (ACh).

Treated with different concentrations of acetylcholine, the cells reliably produced currents in the nanoampere range (Fig. 2).

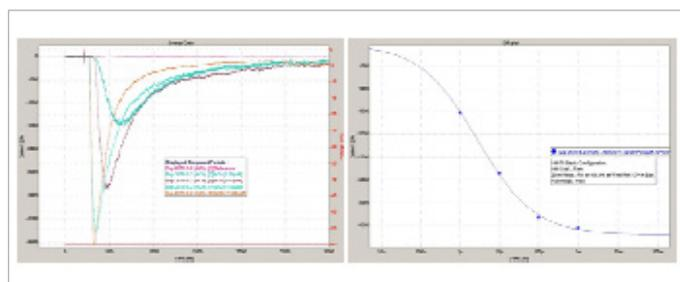
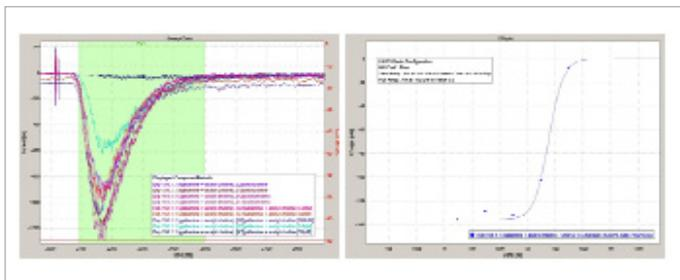


Fig. 2: Left: raw traces with increasing ACh concentration. Right: Agonist (ACh) Hill fit.

Exposure to an increased concentration of ACh showed dose dependency with an  $EC_{50}$  of  $2.67 \pm 1.1 \mu M$  ( $n=11$ ) ( $7.8 \mu M$ , 2) (Fig. 2).

Using cells directly from the freezer gives a lot of flexibility and less day to day variation. In this set of experiments, the frozen cells were used in an antagonist concentration-response experiment using gallamine, a blocker of nAChR $\alpha 1$  (3), to investigate whether the obtained current was directly mediated via the acetylcholine receptor  $\alpha 1$ .

Gallamine blocked the current in a concentration-dependent manner and had an  $IC_{50}$  value of  $5.32 \pm 0.25 \mu M$  ( $n=5$ ) (Fig. 3).



**Fig. 3:** Left: Raw current traces with increasing concentrations of gallamine with constant ACh concentration. Right: Hill fit on gallamine data.

## Methods

The cells are grown according to the Sophion standard operating procedure (SOP) for RD(TE671) cells. Physiological Ringer's solution was used on QPatch. When washing, extracellular Ringer's solution with 2 unit/ml acetylcholinesterase was used for washing.

## Assay

A whole cell protocol modified from a standard CHO protocol was used to obtain the whole cell configuration. During the addition of acetylcholine, the cells were kept at two different potentials, -60 and -90 mV and the channels were opened with a 200-sec interval.

## Author:

Rasmus B Jacobsen, Application scientist

Sophion Bioscience A/S, Baltorpevej 154, 2750 Ballerup, Denmark  
Phone: +45 4460 8800, E-mail: info@sophion.com

sophion.com

## Conclusion

These experiments demonstrate that the QPatch is an ideal platform for reliable dose-response experiments with a fast ligand-gated channel. Using a specific agonist and antagonist, it was confirmed that the observed currents indeed are due to the activation of acetylcholine receptors.

The workload involved in creating a cell line expressing a specific gene of interest can sometimes set the limit for initiating a set of pilot experiments. To be able to work with receptors being endogenously expressed in well-known cell lines can be not only cost-reducing but also instrumental in running initial studies of certain receptor targets. We have in this study examined the RD(TE671) cell line for its endogenous expression of nAChR $\alpha$ 1 and validated its capability to be used on the QPatch platform. We conclude that the overall success rates in obtaining stable seals and completed experiments are at a level that clearly identifies the RD(TE671) assay as feasible and meaningful on the QPatch.

## References:

1. Morris A et al. 1991. Two isoforms of the muscle acetylcholine receptor alpha-subunit are translated in the human cell line TE671. *FEBS Lett* 295:116-8
2. Shao Z et al. 1998. Potentiation and inhibition of nicotinic acetylcholine receptors by spermine in the TE671 human muscle cell line. *J Pharmacol Exp Ther* 286:1269-76
3. Benson JA. 1992. Electrophysiological Pharmacology of the Nicotinic and Muscarinic Cholinergic Responses of Isolated Neuronal Somata from Locust Thoracic Ganglia. *J Exp Biol* 170: 203-233