# PARALLEL PATCH CLAMP OF $\alpha$ 7 NICOTINIC ACETYLCHOLINE RECEPTOR CHANNELS

<sup>1</sup>SØREN FRIIS

<sup>1</sup>Sophion Bioscience A/S, 154 Baltorpvej, DK-2750 Ballerup, Denmark.

The human  $\alpha$ 7 nicotinic acetylcholine receptor  $(\alpha$ 7-nAChR) is a neuronal ligand-gated, fast desensitizing, non-selective cation channel. It is pentahomomeric, consisting of five 50 kD  $\alpha$ 7 subunits, each composed of 502 amino acids. The  $\alpha$ 7-nAChR is involved in memory and cognition, and it is widely study addresses the possibility of efficient and distributed throughout the nervous system, especially in cholinergic neurons projecting to hippocampus and cortex. The  $\alpha$ 7-nAChR has also been found to be associated with pathophysiological states. Importantly, it is involved in widespread human neuro-degenerative and psychotic disorders, including Alzheimer's disease and schizophrenia. Therefore the therapeutic potential of  $\alpha$ 7-nAChR

## MATERIALS AND METHODS

**Cells:** For the studies were used GH4-C1 cells expressing rat  $\alpha$ 7-nAChR. The cells were growth-arrested 48 hours prior to use with 0.5 mM Na-Butyrate, detached from the flask with trypsin and sub-sequently kept in culture medium according to Sophion standard procedures. The isolated cells in solution appeared spherical. For a large number of these experiments frozen vials of cells were thawed just before experiment execution and used directly on the QPatch. This has been found to be an ideal to way to maintain a stable supply of suitable cells for this and other cell lines.

Ringer's solutions: Extracellular (mM): 145 NaCl, 4 KCl, 10 HEPES, 2 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>. Intracellular (mM): 120 KF, 20 KCl, 0.1 EGTA, 10 HEPES,  $2 \text{ MgCl}_2$ .



is substantial, and electrophysiological and pharmacological characterization of the receptor has become an increasingly important issue. The fast kinetics (milliseconds) makes precise patch clamp measurements a challenging task. The present precise characterization of  $\alpha$ 7-nAChR function and pharmacology by means of the QPatch automated patch clamp system. Compared to conventional whole-cell patch clamp the system enables a highly increased throughput by simultaneous and asynchronous operation of 16 parallel patch clamp experiments.

**Compound application:** We used the QPlates XL that hold 250  $\mu$ L in the waste. This allowed execution of advanced ligand- and compound application protocols including multiple applications. Acetyl choline (ACh) and nicotine were used as agonists. Agonist and test compound application protocols were defined in the QPatch assay software. Eight teflon coated pipettes applied the compounds via pipetting wells and the QPlate integrated glass flow channels. During recording, the membrane potential was held at -90 mV.

Sealing rates: The whole-cell success rate was about 80%. However, the final success rate with useful completed experiments was less due to lack of expression in all cells.

## Figure 1. Current rise time

Addition of ligands, either ACh or nicotine, elicited typical lpha7-nAChR currents with rise times of 5-10 ms (5 - 95 %). Here the current is induce by 10 mM ACh. The rise time reflects the solution exchange rate rather than the actual opening rate of the  $\alpha$ 7-nAChR.



## Figure 2. ACh esterase eliminates residual ACh

Multiple additions of the same ligand concentration ( 1 mM ACh) led to decreasing amplitudes of peak lpha7-nAChR currents, as seen from the 4 sequential raw current recordings (A) and the resulting I-t plot (B). Because the concentration-response graph for desensitization is left-shifted to that of activation, even minute agonist concentrations remaining after washes will partially desensitize  $\alpha$ 7-nAChR. Enzymatic removal of the ACh with ACh esterase reduced the problem, as seen from the I-t plot presenting 8 successive applications of 1 mM ACh (**C**). Green points represent leak currents.



Between each application is a 150 sec resting period.

## <sup>1</sup>MORTEN R. SUNESEN

<sup>2</sup>Sophion Bioscience, Inc., North Brunswick, NJ 08902, USA

<sup>2</sup>CHRIS MATHES



## Figure 5. Assay reproducibility and stability

1000 1500 2000 2500 3000 Time Isec1

A: Two ACh concentration-response experiments were performed with a time interval of approximately 45 minutes. Green points represent leak currents. **B**: The two Hill fits were practically identical demonstrating a high degree of reproducibility. **Table**: Comparison between mean peak  $\alpha$ 7-nAChR current amplitudes of 5 experiments similar to the one presented in the panel A. At no ACh concentration is the ratio between the two time-separated peak current amplitudes significantly different from 1.00.

1E-1 1E0 1E1 1E2 1E3 1E4 1E5 1E6 Concentration [μΜ]



## Figure 6. Effect of Nicotine - partial agonism

Concentration-response analysis for the effect of nicotine on  $\alpha$ 7-nAChR peak currents. Trace **A** shows the response to 300  $\mu$ M ACh for comparison. Eight increasing concentrations (traces **B-I**) of nicotine were employed.



A and C: AUC-t plot and Hill fit for the effect of increasing concentrations of nicotine on AUC. B and D: I-t plot and Hill fit for the effect of nicotine on peak  $\alpha$ 7-nAChR currents. EC<sub>50</sub> value compare well to values obtained with Dynaflow (EC<sub>50</sub>=9.9  $\mu$ M).



### <sup>3</sup>JOHN DUNLOP <sup>3</sup>MARK BOWLBY

<sup>3</sup>Discovery Neuroscience, Wyeth Research, CN-8000, Princeton, NJ 08543, USA.

Figure 4. ACh: Peak vs. area under curve (AUC) analysis A: Concentration-response relationship for the effect of ACh on  $\alpha$ 7-nAChR current amplitudes. The graph led to EC<sub>50</sub> = 2.0±0.7 mM (n=8). Similar recordings with the ultra-fast solution exchange system Dynaflow led to  $IC_{50} = 0.38$  mM (data kindly provided by Wyeth Research) indicating that the 5-fold increased QPatch IC<sub>50</sub> value is a result of a slower solution exchange.

**B**: An alternative 'area-under-curve' (AUC, i.e. the integrated charge transfer) analysis led to  $IC_{50} = 26 \pm 7 \mu M$  (n=6) which compares well to the value obtained with Dynaflow, 29  $\mu$ M.

[ACh]	(peak 1/peak 2)·100 (% ± SEM)
41 µM	101 ± 13
123 µM	105 ± 8
370 µM	107 ± 10
1.11 mM	101 ± 11
3.33 mM	87 ± 12





agonist (300 µM ACh) response (control). **B**: Original current recording of NS1738-stimulated  $\alpha$ 7-nAChR current. The NS1738 concentration was  $30 \,\mu$ M.

## CONCLUSION

Despite fast activation and subsequent desensitization which leaves a very brief time window for recordings, it was possible to efficiently study the effects of agonists (full or partial), antagonists and modulators on α7-nAChR expressed in cultured GH4-C1 cells. Current peaks with rise times of 5-10 ms were routinely recorded. The results demonstrate that  $EC_{50}$  and  $IC_{50}$  values for peak currents or charge (AUC) can be measured with accuracy in a highly reproducible manner. Optimization of experimental conditions and unmanned operation makes it possible to obtain higher throughput compared to manual patch clamp.

## <sup>4</sup>NIELS J. WILLUMSEN

