

Application Report

CHO-hERG DUO tested on QPatch in multi-hole mode

An introduction to methodologies that are used when running multi-hole experiments on the QPatch

Summary

This report evaluates the CHO-hERG DUO cell line performance on QPatch when the multi-hole technology was used. It has a specific focus on success rates, stability, pharmacology and biophysical properties of the assay but it also gives an introduction to some of the methodology used when running multi-holes.

Introduction

Formation of a tight seal between cell and chip is crucial in patch clamping. It is, therefore, important to optimize the whole-cell protocol for each cell line of interest. We have found that the use of so-called timed protocols where suction is changed at specific times is a very efficient way to form seals with multi-hole plates.

In multi-hole mode, there are ten cells, which are patched in parallel and there is, therefore, ten times more cell membrane

to rupture when going into whole-cell configuration. We have found that the suction used for whole-cell break-in in multi-hole mode must be harder than in single-hole mode, but still soft enough so that the seal is kept. In Figure 1, two timeline plots are shown and by combining the two plots the total capacitance ($C_{chip}+C_{cell}$) increases 35 pF when performing the whole-cell suction.

Even though the seals are high, one has to take the increased leak into account when using the multi-hole plates. This originates from the fact that ten cells in parallel appear leakier than a single cell because all leak conductance's from the ten holes are summarized.

The QPatch software has several methods for subtracting the ohmic linear leak current. The classic P/n leak subtraction is a method where the cells are exposed to miniature versions of the voltage protocol of use. These miniature sweeps are then used to determinate the leak component, which can then be subtracted from the current. Since the time used for each miniature, the sweep is the same for each voltage protocol, the P/n leak correction is not so practical for assays with long voltage protocols, such as the ones used for hERG measurements, as the total experimental time will potentially affect overall success rates. Other methods must be taken into use.

Online P/n leak subtraction can be seen in Figure 2. The benefit of using this feature is that the leak current is determined between each voltage protocol and this method is, therefore, less sensitive to changes in cell parameters during experiments.

The downside of using this feature with long hERG voltage protocols is the sensitivity for bad sweeps. If the parameters "jump" during one leak sweep, then the whole measured sweep is affected.

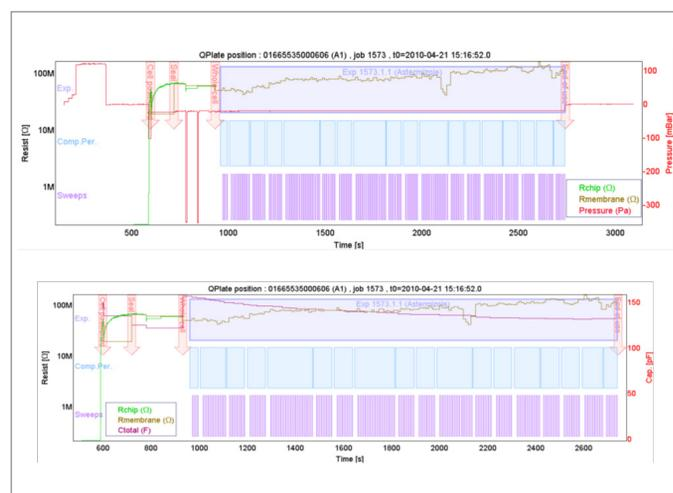


Fig. 1: Timelines from a representative well. By combining the two graphs, the total capacitance increases (red curve bottom plot) when whole-cell suction is applied (red curve top plot).

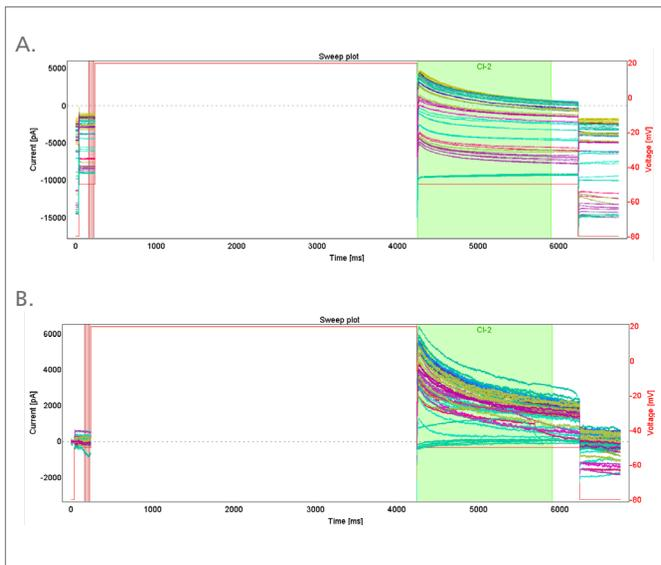


Fig. 2: Example of sweeps from a multi-hole experiment without (a) and with (b) online P/n leak subtraction.

Sweep subtraction is a faster leak method where a specific control antagonist is applied to the cell at the end of the experiments and the current which is left is then subtracted from all other sweeps. The disadvantages of this method are that the seal/leak must stay constant for the whole experiment duration and that the control antagonist must be specific for the current of interest.

Yet another leak subtraction method is available in the QPatch software packages. This is a pre-pulse leak subtraction method where a small pre-pulse or ramp is used for an AC measurement of the leak current at two given potentials thereby allowing for a more accurate leak subtraction at other potentials. This method is therefore mainly useful for IV protocols (Figure 3).

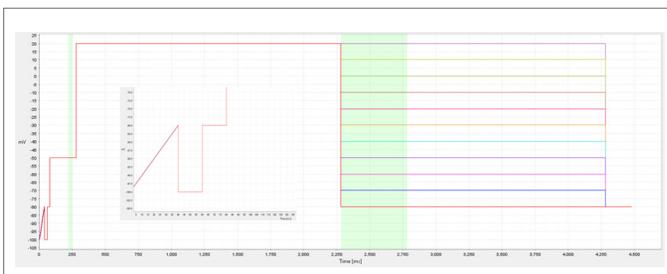


Fig. 3: Voltage protocol with ramp and prepulse for reversal potential investigations with leak corrections.

All three methods available in the software have been tried in this study and they all have their strengths.

In Figure 4 an example of a single experiment/measurement site can be seen, with sweep subtraction and without sweep subtraction. Note that the baseline current is shifted to 0 pA after sweep subtraction.

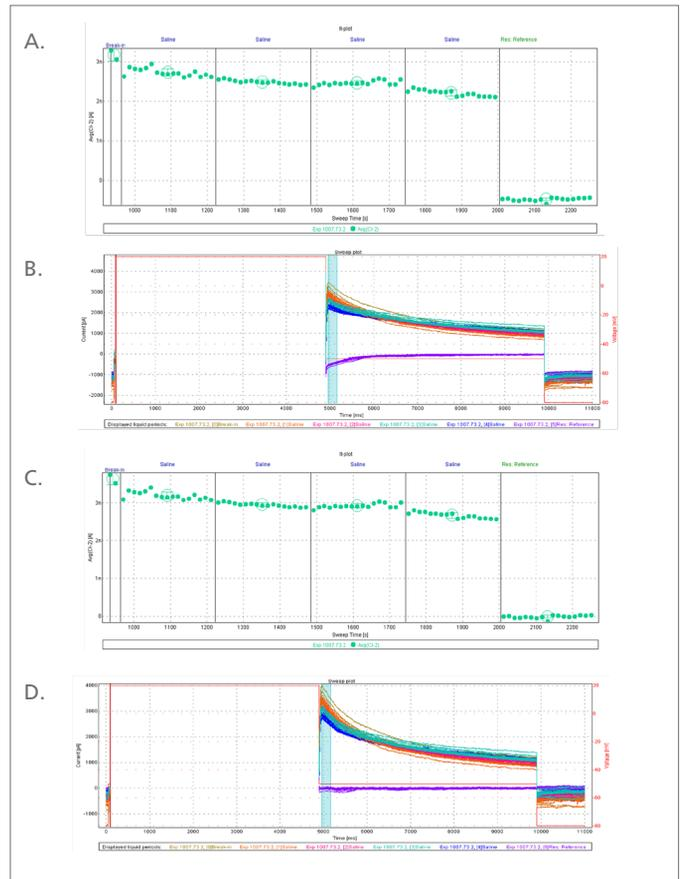


Fig. 4: Example of use of sweeps subtraction, all four figures originate from the same raw data. a and c) current versus time plot (I/t-plot) with and without sweep subtraction respectively b and d) hERG responses with and without sweep subtraction.

Results and discussion

Throughput

The throughput was evaluated in a series of 6 pt. dose-response experiments. Throughput was calculated as the number of experiments that were found useful after analysis i.e. the rightmost column in the histogram (Figure 5).

IC₅₀ values were determined by the use of group Hill fit and each useful liquid period was included in the analysis in the multi-hole experiments (Figure 5).

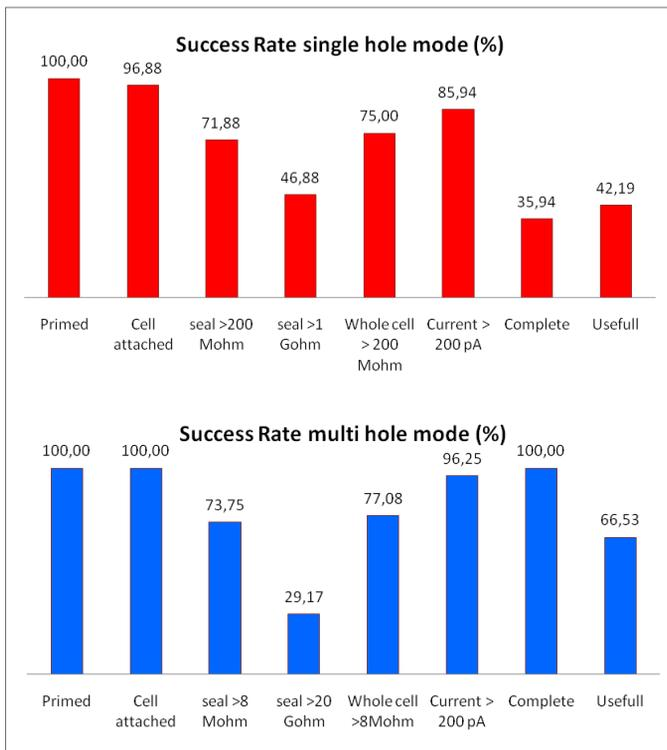


Fig. 5: Top: Success rate in single hole mode (n=64). Bottom: Success rate in multi-hole mode (n=240).

Stability of recordings

Rundown can be a problem when running hERG. The effect on the rundown and current stability was evaluated by measuring the coefficient of variance (CV) and rundown of the last saline period before the compound was added.

CV was measured as $SD/average$ current derived with the assay software from 500 sites on both single-hole and multi-hole QPlates.

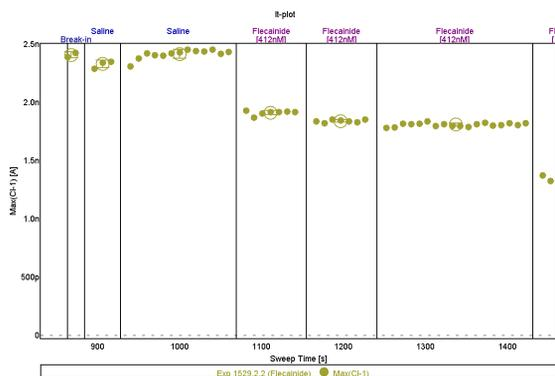


Fig. 6: hERG statistical analysis It-plot. SD and average of current in the last saline period are used for CV determination.

The distribution of the found CV's is shown in Figure 7. CV was found to be slightly but not significantly lower with multi-hole compared to single-hole mode.

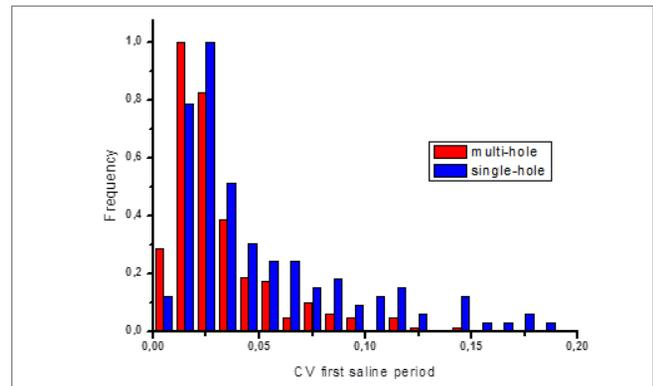


Fig. 7: Distribution of saline CV in single-hole and multi-hole.

The rundown in these hERG experiments could be described with a linear function. The assay software was therefore used to make a linear fit to the last saline period and the rundown rate (% rundown/min) was derived from this result. The distributions of rundown rates are shown in Figure 8. It was found that the rundown was reduced slightly in multi-hole mode, however, it was also found that there was more run-up in multi-hole mode than seen in single-hole mode.

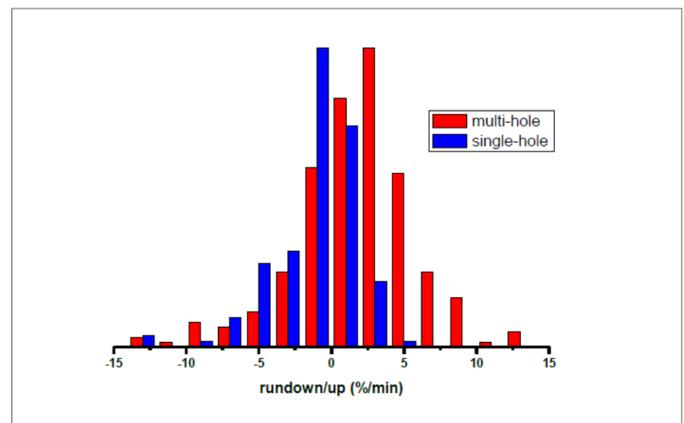


Fig. 8: Distribution of rundown rates in multi-hole and single-hole mode.

Biophysics of hERG in multi-hole mode

When performing current vs. potential experiments (IV-experiments) on reversal potential, an off-line leak protocol is needed. By using a pre-pulse or a pre-ramp, it is possible in the QPatch assay software to subtract a leak at any potential (Figure 3). IV curves by using this method can be seen in Figure 9 when reversal potential is measured to -80 mV.

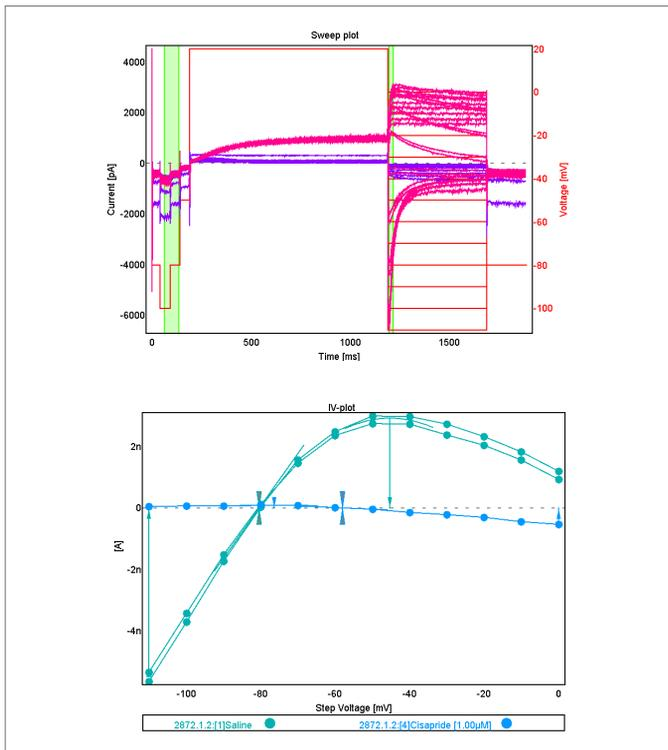


Fig. 9: The current-voltage relationship for determination of the reversal potential. Left: voltage protocol and raw current traces. Right: IV relationship, green line with physiological Ringer's solutions and blue line in the presence of 1 μ M Cisapride.

Pharmacology

We have tested six known blockers of the hERG current. The obtained results were used to validate the multi-hole technology against single-hole recordings and literature values.

Table 1: Half maximal inhibitory concentration (IC_{50}) values from six known blockers of the hERG current. All values are in nM. *) determined by individual Hill fits #) data normalized and IC_{50} value determined by group Hill fit.

	Literature values (nM)	Single-hole mode* (nM)	Multi-hole mode#
Astemizole	37 ⁽²⁾	23 (n=8)	16 (n=153 data points)
Cisapride	20-10,000 ⁽³⁾	56 (n=9)	26 (148 data points)
Flecainide	3,910 ⁽⁴⁾	1360 (n=1)	1779 (97 data points)
Pimozide	2-20 ⁽³⁾	41 (n=2)	13 (103 data points)
Quinidine	200-2,000 ⁽³⁾	1259 (n=4)	945 (108 data points)
Verapamil	100-1,000 ⁽³⁾	340 (n=3)	137 (107 data points)

Conclusion

By using an appropriate whole-cell protocol and a proper leak subtraction the multi-hole technology is a powerful tool in testing hERG pharmacology.

We have been able to increase the throughput due to more experiments being accepted both because more experiments complete (42 vs. 66 %, Figure 5), but also because the stability of the recording is improved (Figure 6).

The pharmacological data are in good agreement with both single-hole experiments and already published literature values.

Methods

CHO-hERG DUO cells were cultured and harvested according to Sophion SOP.

References

1. Armstrong, CM, Bezanilla, F. Charge movement associated with the opening and closing of the activation gates of the Na channels. *J.Gen. Physiol.* 1974. 63:533-552.
2. Brown, AM. hERG block, QT liability and sudden cardiac death. *Novartis.Found.Symp.* 2005. 266:118-131.
3. Ficker, E, Dennis, A, Kuryshv, Y, Wible, BA, Brown, AM. hERG channel trafficking. *Novartis.Found.Symp.* 2005. 266:57-69.
4. Kuryshv, YA, Ficker, E, Wang, L, Hawryluk, P, Dennis, AT, Wible, BA, Brown, AM, Kang, J, Chen, XL, Sawamura, K, Reynolds, W, Rampe, D. Pentamidine-induced long QT syndrome and block of hERG trafficking. *J.Pharmacol.Exp.Ther.* 2005. 312:316-323.