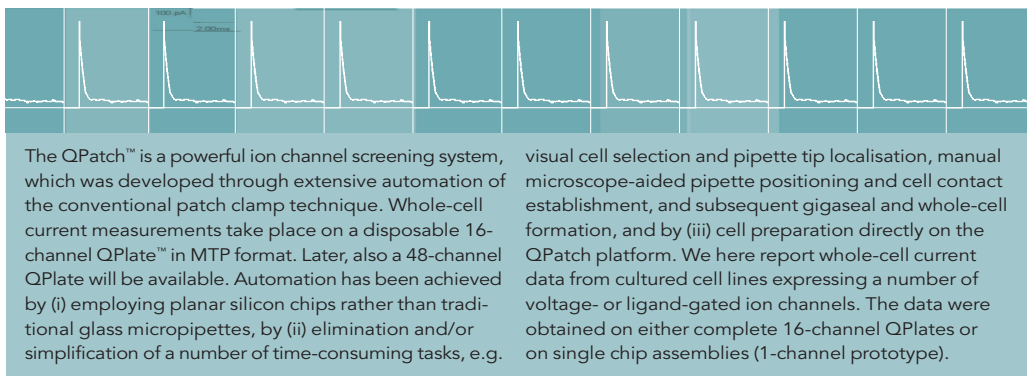


# AUTOMATION OF THE PATCH-CLAMP TECHNIQUE: TECHNICAL VALIDATION THROUGH CHARACTERIZATION OF VOLTAGE-GATED POTASSIUM CHANNELS AND LIGAND-GATED ION CHANNELS

KAREN KRZYKOWSKI | RIKKE L. SCHRÖDER | TRINE LJUNGSTRØM | JONATAN KUTCHINSKY | SØREN FRIIS | MARGIT ASMILD | RAFAEL TABORSKI | SIMON PEDERSEN | RAS K. VESTERGAARD | RASMUS B. JACOBSEN | NATHALIE HÉLIX | CLAUS B. SØRENSEN | MORTEN BECH | JEPPE K. CHRISTENSEN | SIMON DUBEAU | NIELS J. WILLUMSEN



**MATERIALS**  
**ION CHANNEL PROTEINS**  
 For patch clamp measurements were used a number of different ion channels: Voltage gated ion channels (VGIC): HERG, KCNQ4. Ligand gated ion channels (LGIC): GABA<sub>A</sub>-R, nACh-R, ASIC

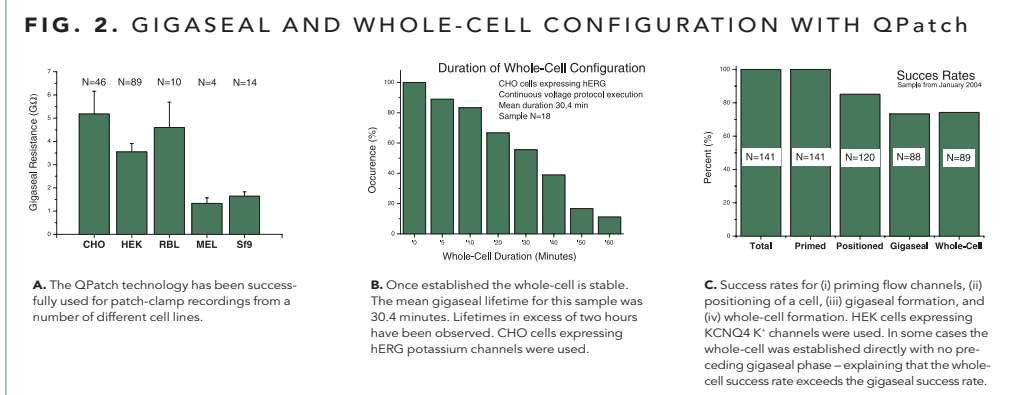
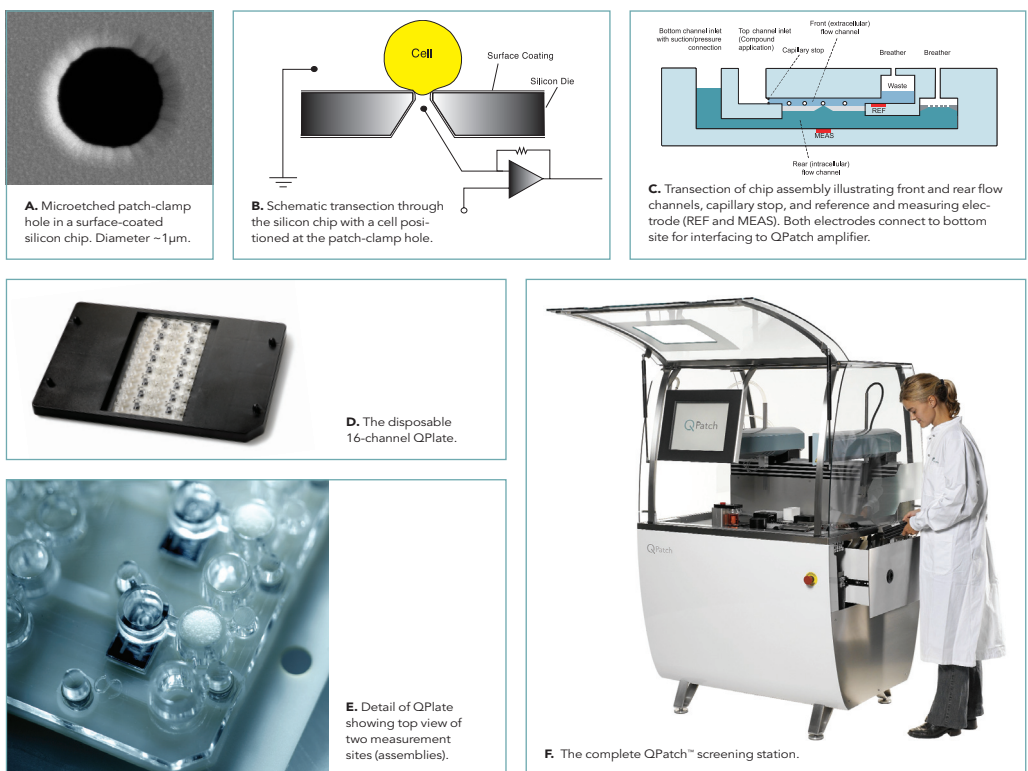
**CELL CULTURES**  
 The ion channels were expressed in either HEK-293 or CHO-k1 cells. These cell lines were grown according to standard culturing protocols (Kutchinsky et al., Assay and Drug Development Technologies 1:685-693, 2003). All cells were grown at 37°C. After harvest, cells could be kept for up to 4 hours in the cell storage facility on the QPatch platform with no change in quality or ability to form gigaseals.

visual cell selection and pipette tip localisation, manual microscope-aided pipette positioning and cell contact establishment, and subsequent gigaseal and whole-cell formation, and by (iii) cell preparation directly on the QPatch platform. We here report whole-cell current data from cultured cell lines expressing a number of voltage- or ligand-gated ion channels. The data were obtained on either complete 16-channel QPlates or on single chip assemblies (1-channel prototype).

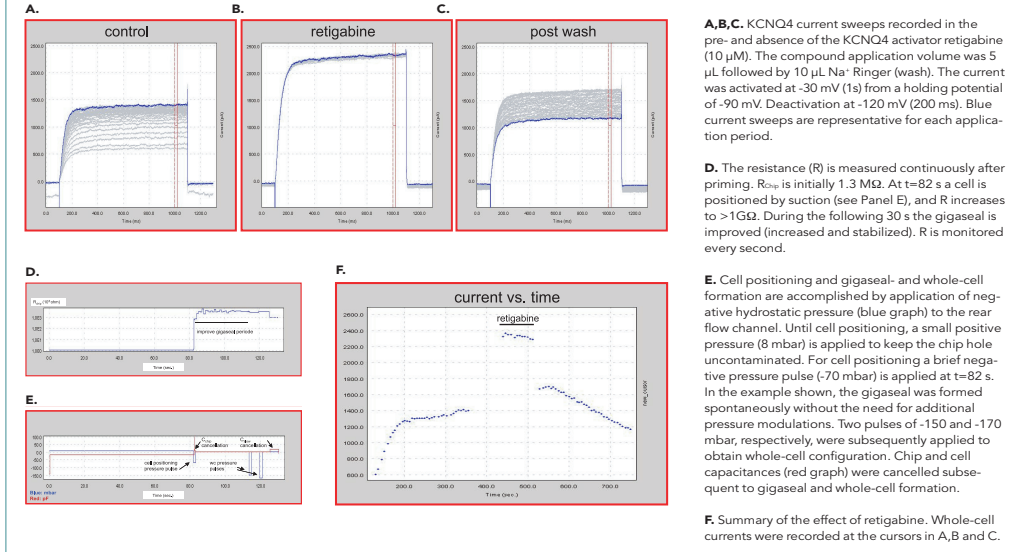
**ELECTROPHYSIOLOGY**  
 Whole-cell measurements were done either on single planar silicon chips or on complete 16-channel QPlates (Figure 1D; Kutchinsky et al., 2003) with either a 16-channel Sophion patch clamp amplifier or with a HEKA (EPC-9, HEKA Elektronik, Germany) commercial amplifier. The chips had micro-etched patch clamp holes with diameters of approximately 1 μm (Figure 1A) and resistances of 2.04±0.02 MΩ (N=274) in symmetrical physiological saline.

**SOLUTIONS AND DRUGS**  
 The physiological Ringer solutions consisted of (in mM). Extracellular Na<sup>+</sup> Ringer: 140 NaCl, 4 KCl, 2 CaCl<sub>2</sub>, 1 MgCl<sub>2</sub>, 10 HEPES (pH 7.4). Intracellular K<sup>+</sup> Ringer: 120 KCl, 5.4 CaCl<sub>2</sub>, 1.8 MgCl<sub>2</sub>, 30/10 KOH/EGTA, 10 HEPES, 4 ATP, 0.4 GTP (pH 7.2). Verapamil was from Sigma, Switzerland. rBeKm-1 was from Alomone Labs, Israel. GABA and retigabine were from NeuroSearch, Denmark. In LGIC experiments solutions were added in 2 μL volumes by an autosampler.

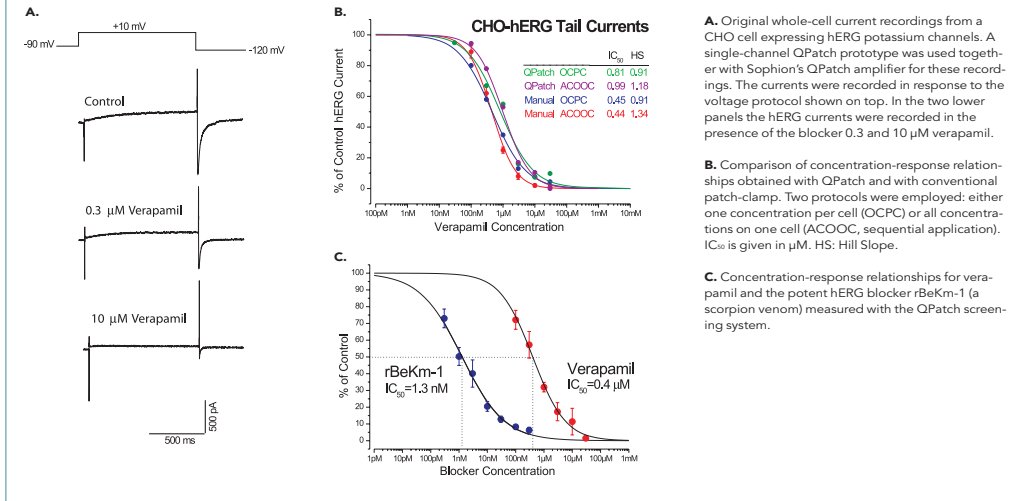
**FIG. 1. THE QPatch PATCH-CLAMP SCREENING SYSTEM**



**FIG. 2. GIGASEAL AND WHOLE-CELL CONFIGURATION WITH QPatch**

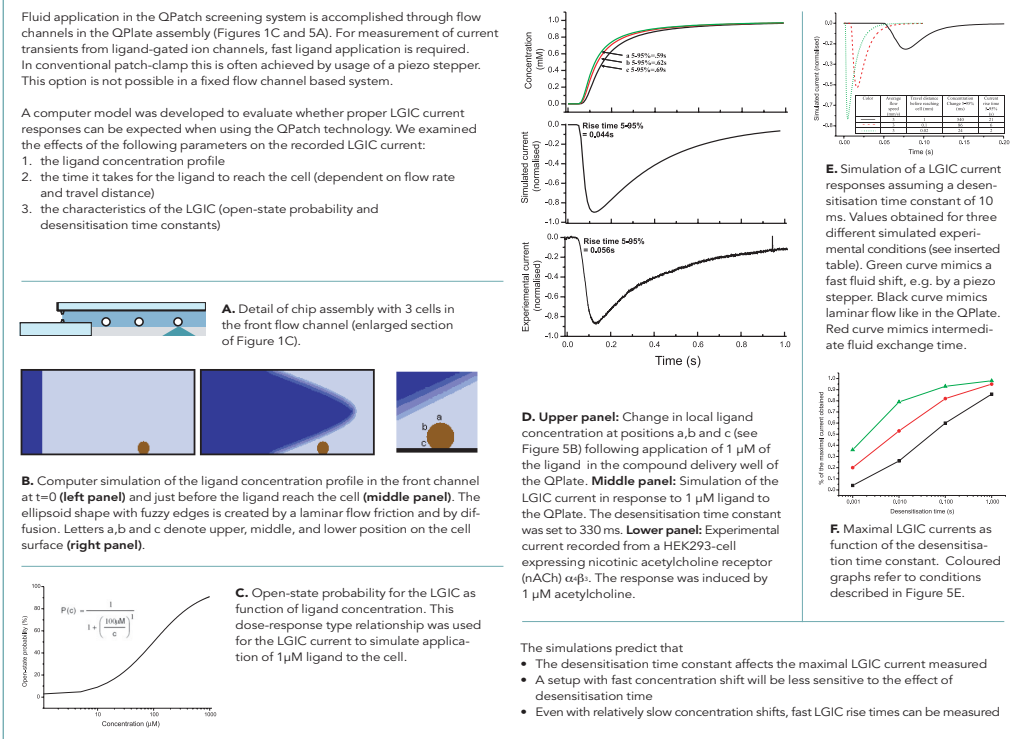


**FIG. 3. SCREENING WITH QPatch: MEASUREMENTS OF KCNQ4 CURRENTS ON A 16-CHANNEL QPlate**

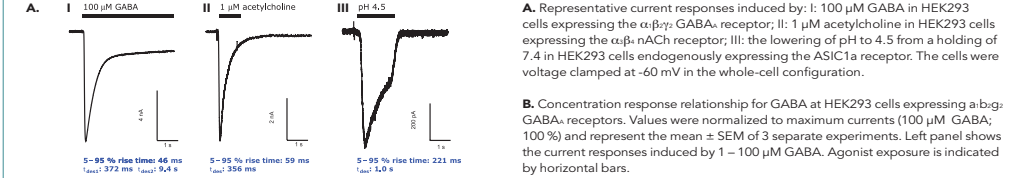


**FIG. 4. PATCH-CLAMPING WITH QPatch: CONCENTRATION-RESPONSE RELATIONSHIPS FOR hERG CHANNEL BLOCKERS**

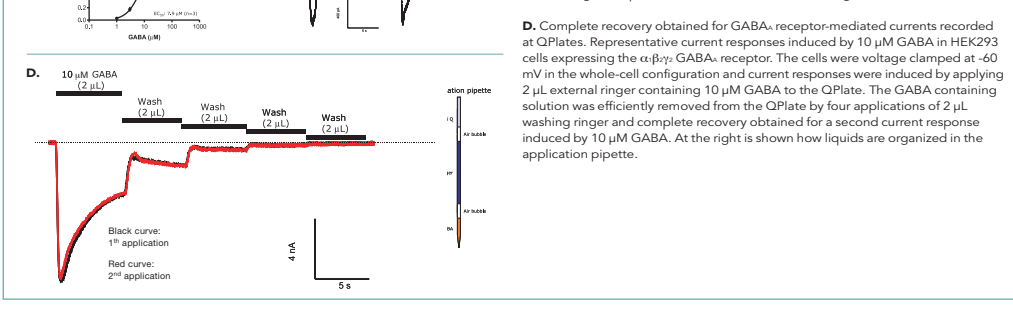
**FIG. 5. CAN QPatch MEASURE LIGAND-GATED ION CHANNEL CURRENTS? COMPUTER MODEL PREDICTIONS**



**FIG. 5. CAN QPatch MEASURE LIGAND-GATED ION CHANNEL CURRENTS? COMPUTER MODEL PREDICTIONS**



**FIG. 6. RECORDING OF LIGAND-GATED ION CHANNEL CURRENTS**



**SUMMARY**  
 High-quality whole-cell recordings from a number of ion channels expressed in cultured cell lines have demonstrated that voltage-gated as well as ligand-gated ion channel proteins can be efficiently targeted by the QPatch screening technology. Automation of the patch-clamp technique will substantially increase throughput in future ion channel drug discovery