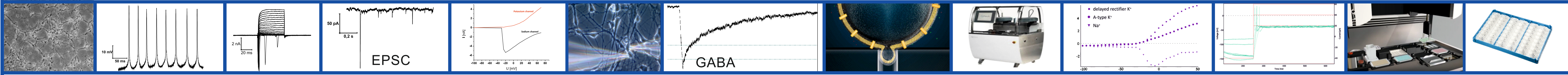


Electrophysiological characterization of human induced pluripotent stem cell-derived dopaminergic neurons using manual and automated patch clamp systems

Denise Franz¹, Jan Gimsa¹, Oliver Klink¹ and Hervør Lykke Olsen²



Introduction

We present a basic electrophysiological characterization of the Dopa.4U, human induced pluripotent stem cell-derived dopaminergic neurons of Axiogenesis (Cologne, Germany). That included examining the voltage-gated ion channels as well as ligand-gated GABA receptors on the manual patch clamp and the automated QPatch system (Sophion - Biolin Scientific, Ballerup, Denmark).

Our results suggest the feasibility of an automated electrophysiological characterization of suspended neuronal cells.

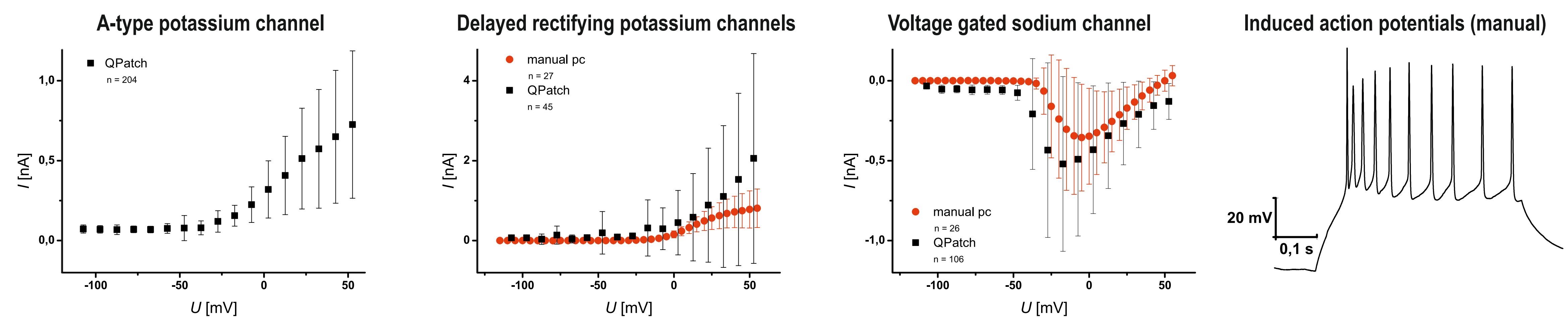
Materials & Methods

The Dopa.4U were cultured on Geltrex[®] or Matrigel[®] coated culture dishes for at least 9 days. Because of using suspended cells for measurements with automated system, the following results consider and compare attempts of detached Dopa.4U. Of three different harvesting-compounds tested, Accutase[®] (Sigma Aldrich) seemed to allow for the mildest treatment. All manual experiments were executed using the triple EPC 10 amplifier (HEKA Elektronik, Germany).

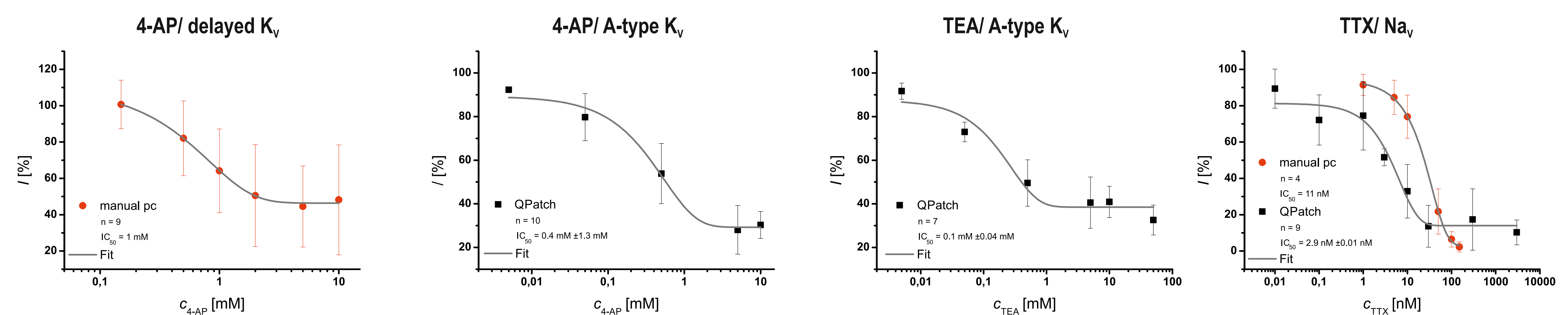
Results

In manual patch clamp experiments, two voltage-gated ion channel types (K_v and Na_v) were characterized with the potassium-channel blocker 4-Aminopyridine (4-AP) and the sodium-channel blocker Tetrodotoxin (TTX) leading to IC_{50} values of 1 mM and 1 nM, respectively. With the automated QPatch system, a gigaseal rate of up to 90 % and a success rate of up to 70 % have been achieved. Detailed analysis of the voltage gated ion channels revealed 83 % cells with rapidly inactivating A-type potassium channels, 43 % cells with sodium channels, 18 % cells with delayed rectifying potassium channels, and 21 % cells with both, A-type and delayed rectifying potassium channels. Pharmacological investigations with different channel blockers provided stable results for the A-type potassium channel with 4-AP ($IC_{50} = 0.4 \text{ mM} \pm 1.3 \text{ mM}$) and Tetraethylammonium (TEA; $IC_{50} = 0.1 \text{ mM} \pm 0.04 \text{ mM}$), as well as for the sodium channel with TTX ($IC_{50} = 2.9 \text{ nM} \pm 0.01 \text{ nM}$). First experiments have shown that the QPatch system allows for reproducing induced action potentials, which were obtained with the manual system in the current clamp mode.

Current-Voltage-Curves



Dose-Response-Curves



Summary

	Gigaseal rate	Success rate	A-type K_v	Delayed K_v	Na_v	GABA currents	Induced AP
manual	≤ 97 %	≤ 96 %	-	$U_{open} = -20 \text{ mV} \pm 8 \text{ mV}$ → 4-AP (with adherent cells) $IC_{50} = 1 \text{ mM}$	$U_{open} = -35 \text{ mV} \pm 8 \text{ mV}$ → TTX (with adherent cells) $IC_{50} = 11 \text{ nM}$	only in adherent cells	possible
automated	≤ 90 %	≤ 70 %	$U_{open} = -32 \text{ mV} \pm 7 \text{ mV}$ → 4-AP $IC_{50} = 0,4 \text{ mM} \pm 1,3 \text{ mM}$ → TEA $IC_{50} = 0,1 \text{ mM} \pm 0,04 \text{ mM}$	$U_{open} = -27 \text{ mV} \pm 14 \text{ mV}$	$U_{open} = -37 \text{ mV} \pm 8 \text{ mV}$ → TTX $IC_{50} = 2,9 \text{ nM} \pm 0,01 \text{ nM}$		possible

Acknowledgment

This research is financially supported by the graduate college *welisa* funded by the DFG. We are very grateful to *Axiogenesis*, especially to Dr. Andreas Ehlich, Dr. Gesa Rascher-Eggstein and Dr. Dietmar Hess, for providing their expertise, and sponsorship in the field of stem cell research. Tusind Tak to *Sophion*, primarily Hervør L. Olsen, Mette T. Christensen, Anders Lindqvist and Naja M. Sørensen, for the opportunity of an internship, support and expert advice in the research area of automated patch clamp systems.