

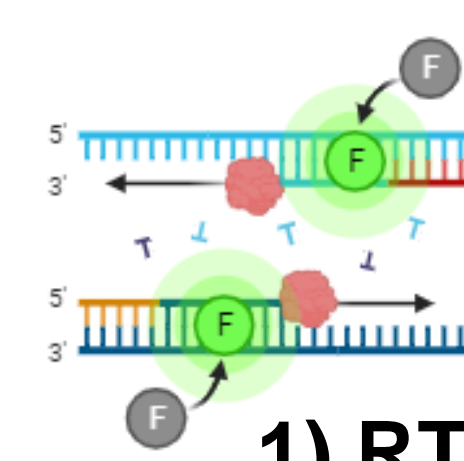
Introduction

K_v7 (KCNQ or M-) channels generate voltage dependent K^+ currents and represent the molecular basis of the M-current (I_M). The five known subtypes ($K_v7.1-7.5$) are characterized by different tissue expression and physiological properties.^[1] In most neurons, native M-channels are composed of $K_v7.2$ and $K_v7.3$ subunits, and are responsible for controlling the neuronal excitability.^[2] During physiological as well as pathophysiological conditions (e.g. epilepsy) channel modulators play a significant role.^[3] Previous studies from our group point out the importance of M-channels in the thalamocortical system, making them interesting targets for further investigation.^[4]

With this project, we thus aim to investigate:

- 1) The developmental expression profile of $K_v7.2$ and $K_v7.3$ in thalamic neurons.
- 2) The modulation of these channels by different compounds including PIP_2 , cholesterol analogues and novel activators.
- 3) The properties of I_M in prethalamic/ thalamic neurons in health and disease.

Methods

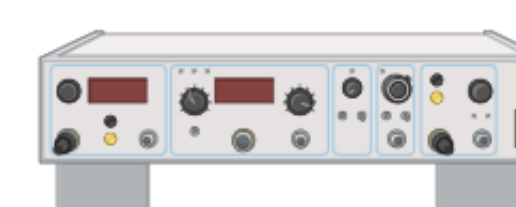


1) RT-qPCR

for the investigation of ion channel expression on mRNA level

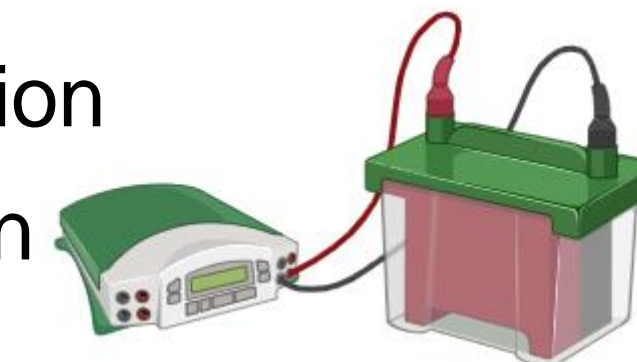
5) Patch-clamp

for the examination of electrophysiological ion channel activity



2) Western blot

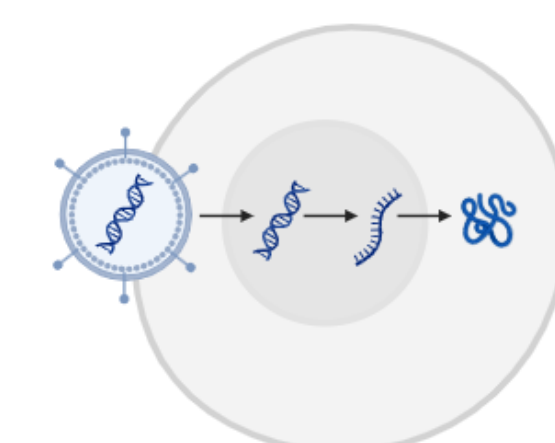
for the investigation of ion channel expression on protein level



Focus on $K_v7.2$ and $K_v7.3$ channels

3) Cell transfection

for generating a cell line with functional M-channel



4) Mouse models

C57BL/6 (wild-type mice)
KCNQ3 KO mice



Figures generated with BioRender

Developmental expression profile of $K_v7.2$ and $K_v7.3$ in mouse brain

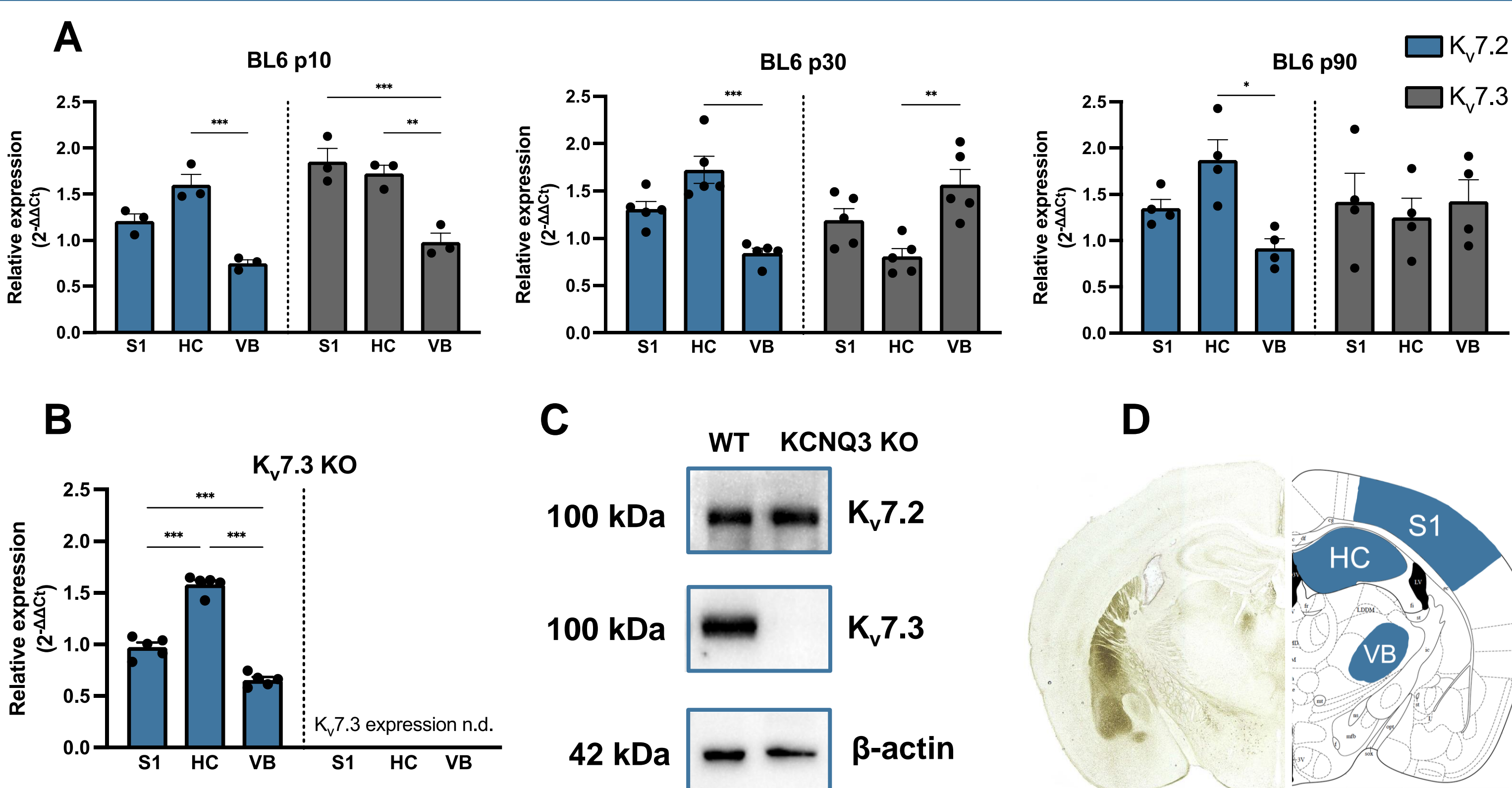


Figure 1: $K_v7.2$ and $K_v7.3$ gene and protein expression in mice during development.

S1: Somatosensory cortex. HC: Hippocampus. VB: Ventrobasal complex. **A)** Relative gene expression in C57BL/6J01aHsd mice. Early infant mice ($n=3$): p10, adolescent mice ($n=4$): p30, adult mice ($n=4$): p90. **B)** Relative gene expression in $KCNQ3$ KO mice ($n=5$): p90 and older. One-way Anova test was applied. Error bars represented as \pm SEM. * $p = 0.033$, ** $p = 0.01$, *** $p < 0.001$. **C)** Representative data for western blotting. $K_v7.2$ and $K_v7.3$ in mouse brain: p30. **D)** Regions of interest in coronal brain slices (Modified from mouse brain atlas, Paxinos and Franklin).

K_v7 modulation with the cholesterol analogue CHIM-L

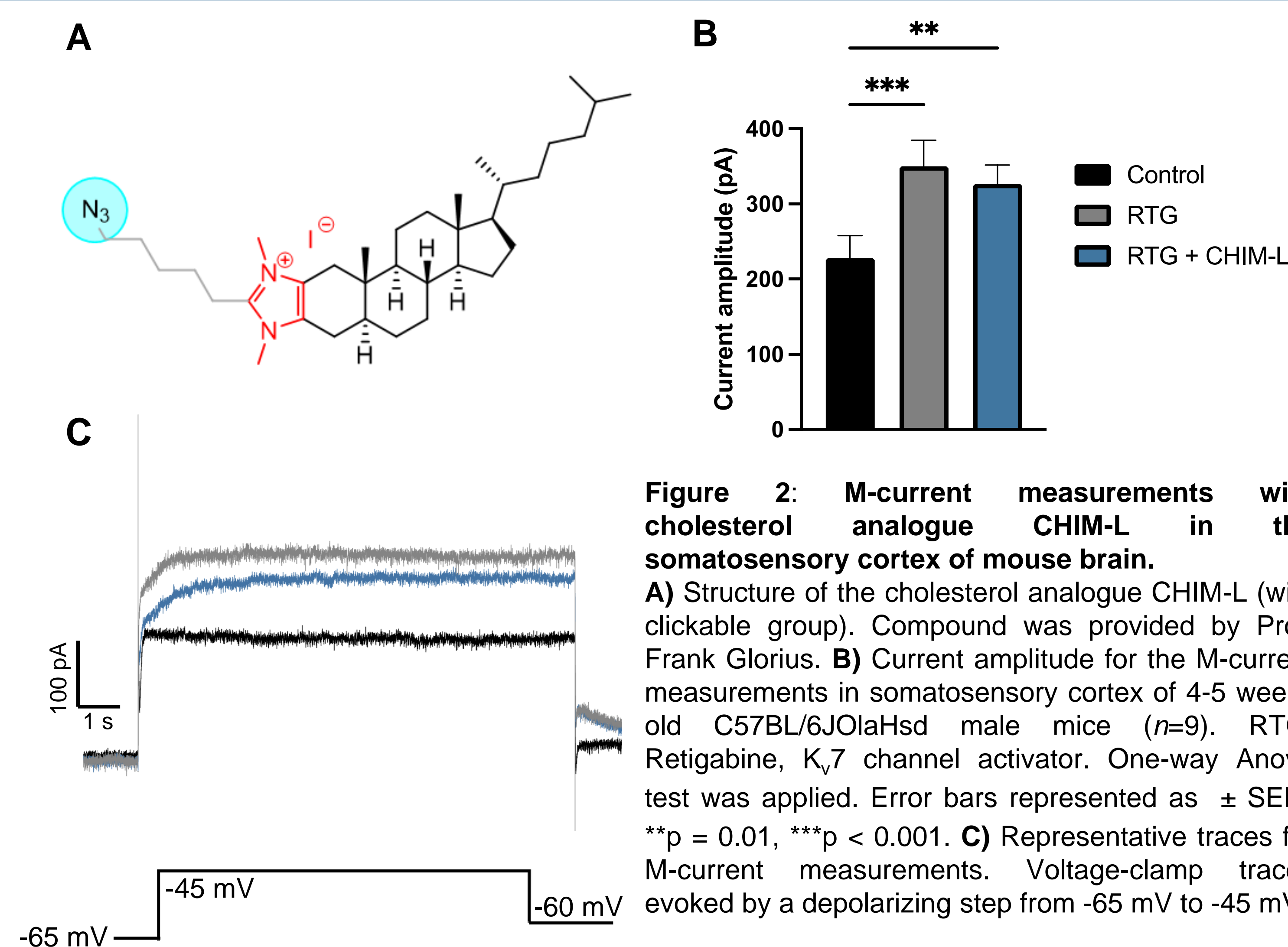


Figure 2: M-current measurements with cholesterol analogue CHIM-L in the somatosensory cortex of mouse brain.

A) Structure of the cholesterol analogue CHIM-L (with clickable group). Compound was provided by Prof. Frank Glorius. **B)** Current amplitude for the M-current measurements in somatosensory cortex of 4-5 weeks old C57BL/6J01aHsd male mice ($n=9$). RTG: Retigabine, K_v7 channel activator. One-way Anova test was applied. Error bars represented as \pm SEM. ** $p = 0.01$, *** $p < 0.001$. **C)** Representative traces for M-current measurements. Voltage-clamp traces evoked by a depolarizing step from -65 mV to -45 mV.

Cell transfection

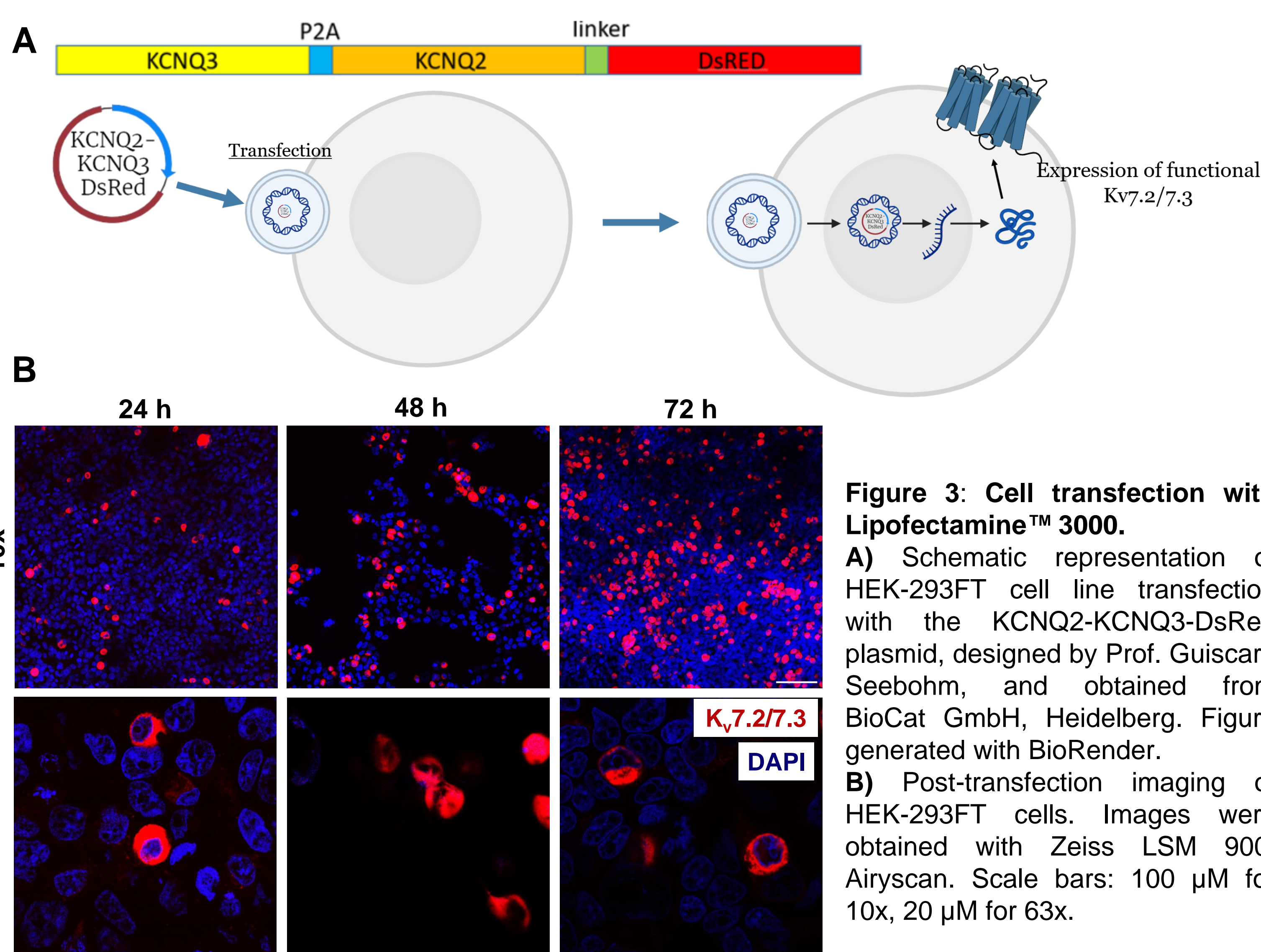


Figure 3: Cell transfection with Lipofectamine™ 3000.

A) Schematic representation of HEK-293FT cell line transfection with the $KCNQ2$ - $KCNQ3$ -DsRed plasmid, designed by Prof. Guiscard Seeböhm, and obtained from BioCat GmbH, Heidelberg. Figure generated with BioRender. **B)** Post-transfection imaging of HEK-293FT cells. Images were obtained with Zeiss LSM 900, Airyscan. Scale bars: $100 \mu\text{m}$ for 10x, $20 \mu\text{m}$ for 63x.

Functionality test of transfected cells

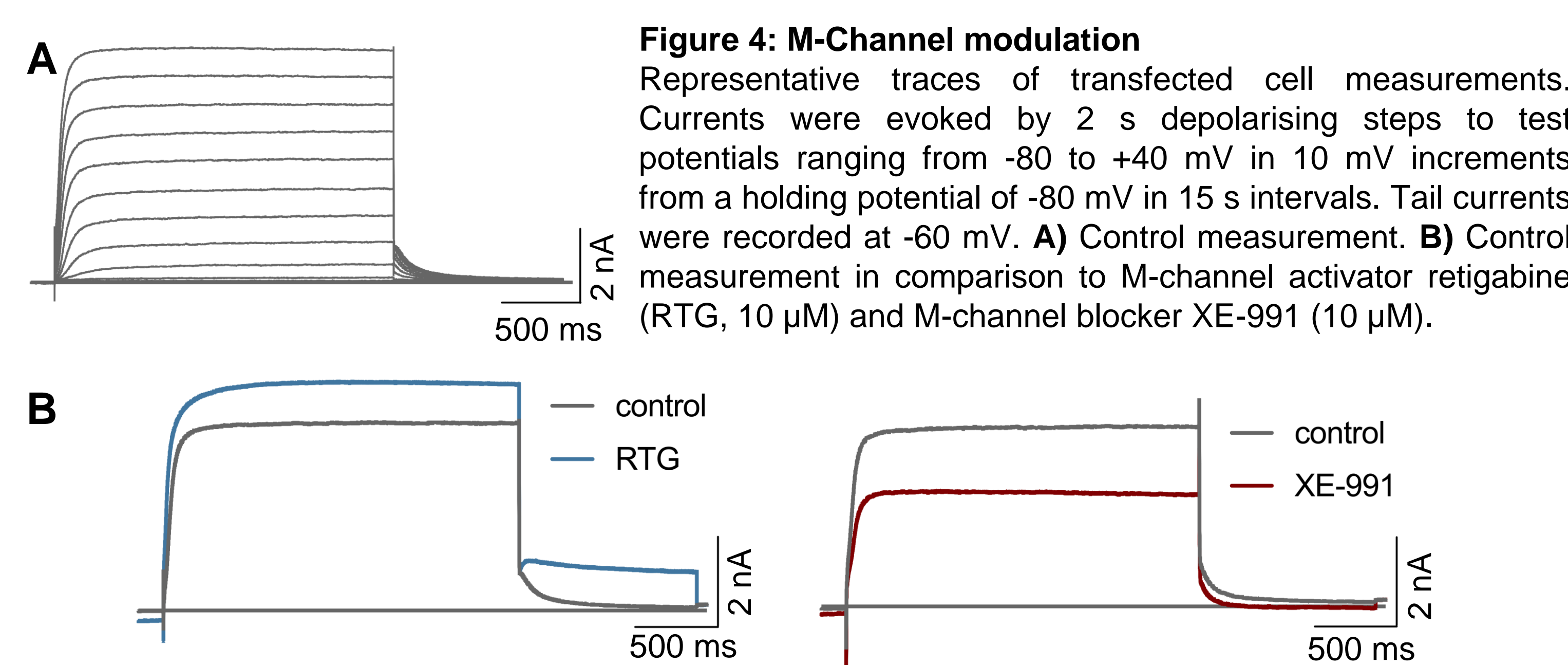


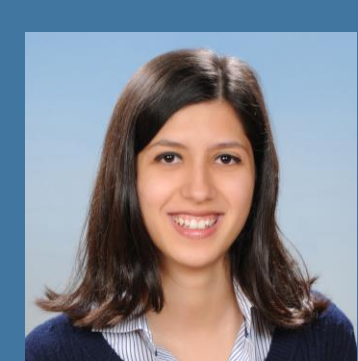
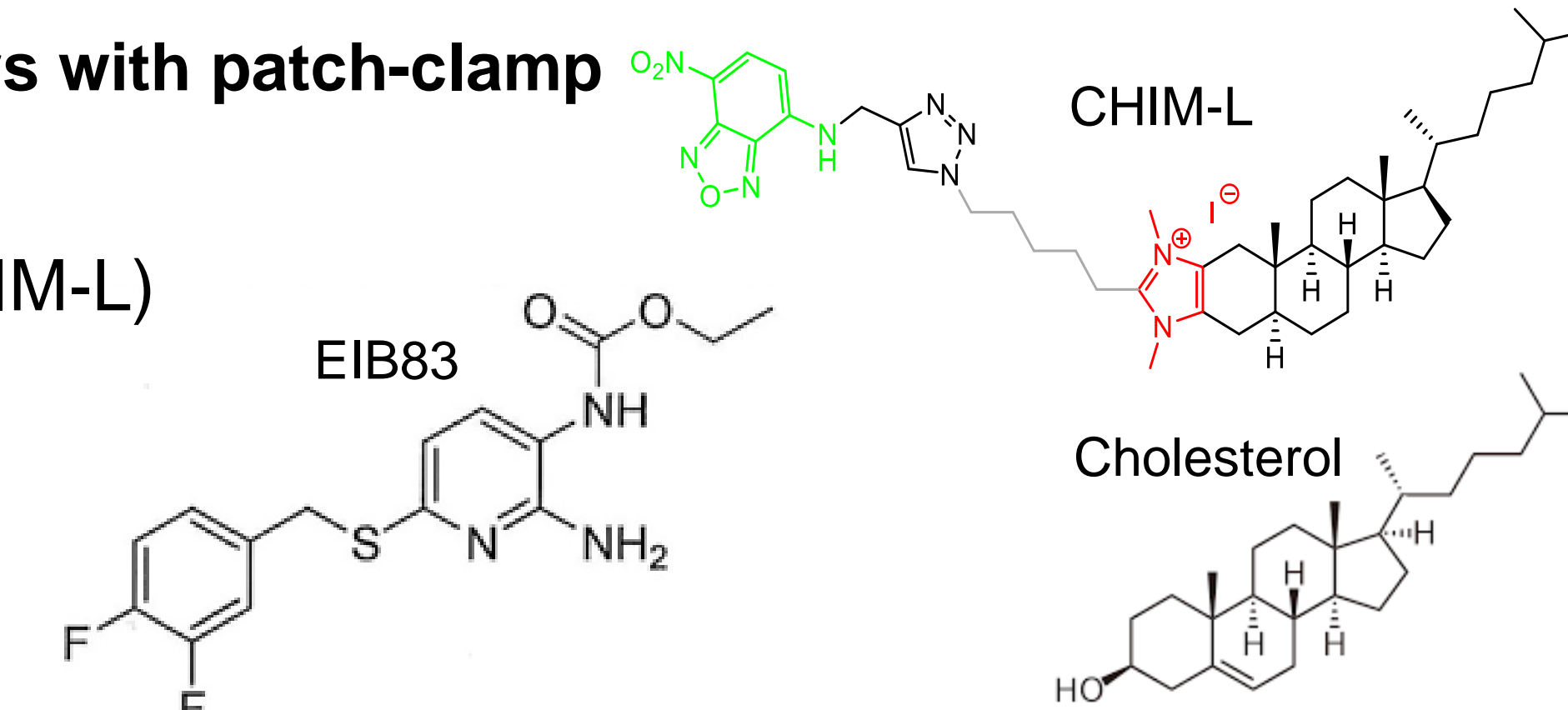
Figure 4: M-Channel modulation

Representative traces of transfected cell measurements. Currents were evoked by 2 s depolarising steps to test potentials ranging from -80 to $+40$ mV in 10 mV increments from a holding potential of -80 mV in 15 s intervals. Tail currents were recorded at -60 mV. **A)** Control measurement. **B)** Control measurement in comparison to M-channel activator retigabine (RTG, $10 \mu\text{M}$) and M-channel blocker XE-991 ($10 \mu\text{M}$).

Future prospects

Test of M-channel modulators with patch-clamp

- Natural cholesterol
- Cholesterol analogues (CHIM-L)
- PIP_2 analogues
- Retigabine analogues



References:

- [1] Dirx *et al.*, *Front. Physiol.*, 2020, S. 1240.
- [2] Brown *et al.*, *Br. J. Pharmacol.*, 2009, 156(8), 1185-1195.
- [3] Borgini *et al.*, *RSC Med. Chem.*, 2021, 12.4, 483-537.
- [4] Cerina *et al.*, *Br. J. Pharmacol.*, 2015, 172.12, 3126-3140.