

Electrophysiological characterization of human dopaminergic neurons derived from LUHMES cells

Udo Kraushaar [1], Dominik Loser [1-4], Timm Danker [2], Clemens Möller [3], Marcel Leist [4]

NMI - Natural and Medical Sciences Institute at the University of Tübingen
Markwiesenstrasse 55
72770 Reutlingen, Germany

E-Mail: udo.kraushaar@nmi.de

[1] NMI Natural and Medical Sciences Institute at the University of Tübingen, Reutlingen, GER; [2] NMI-TT GmbH, Reutlingen, GER; [3] Life Sciences Faculty, Albstadt-Sigmaringen University, Sigmaringen, GER; [4] Doerenkamp-Zbinden Chair for in vitro Toxicology and Biomedicine, University of Konstanz, Konstanz, GER

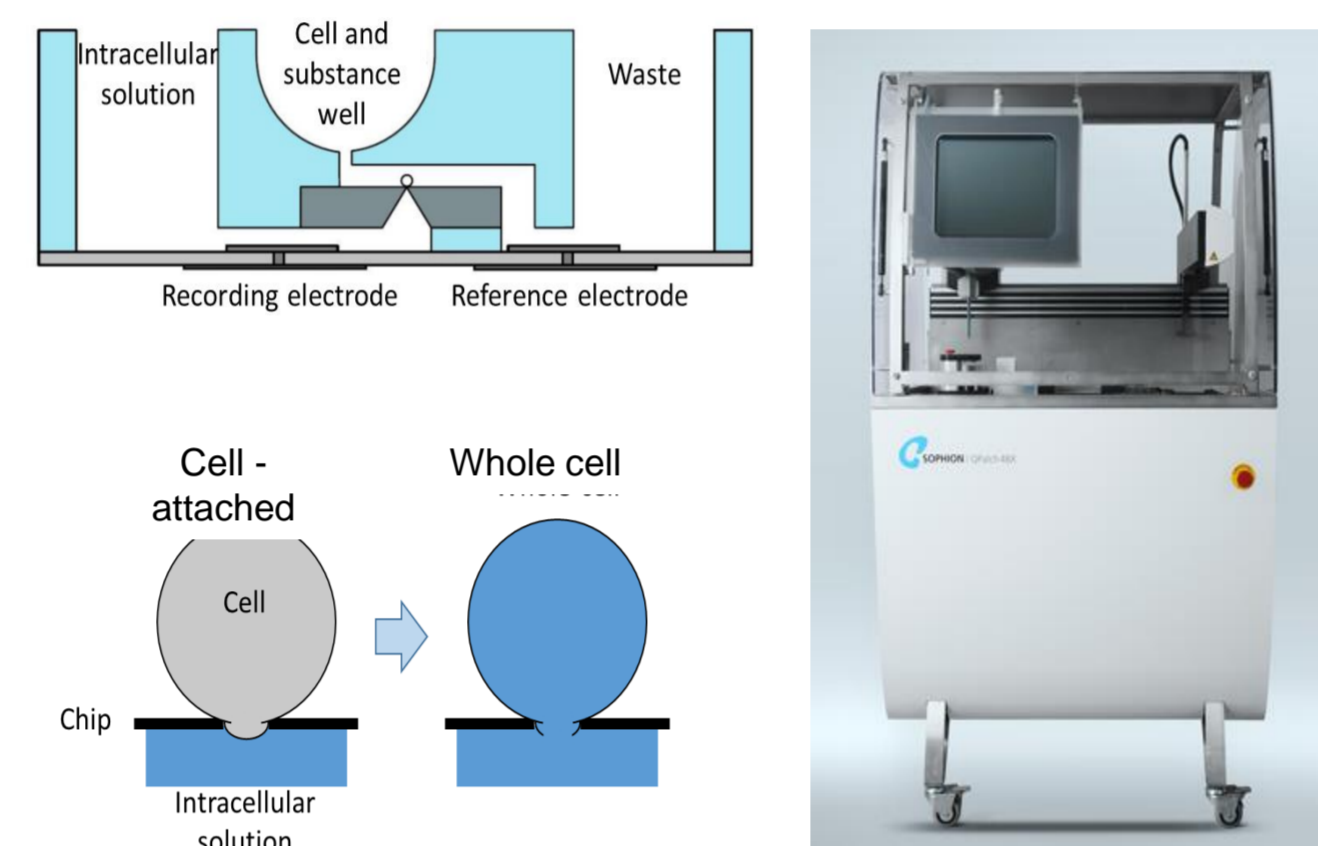
Introduction

The loss of dopaminergic neurons in the substantia nigra plays an important role in the development of the Parkinson's disease. The symptoms of this disease typically occur after around 80 % of these neurons degenerated. This cell decay can be caused or promoted by genetic defects or environmental factors including chemical compounds like pesticides. For a proper testing of neurotoxic effects on these neurons as well as for the development of neuroprotective drugs, assays based on animal primary cells lack pre-

dictivity due to mostly weak correlation between animal and human data. Therefore, models based on human neuronal cells have a high potential to overcome the limitations of animal models. One interesting neuronal cell line is the LUHMES (Lund human mesencephalic) line, which consists of immortalized fetal human mesencephalic cells that can be differentiated into fully post-mitotic dopaminergic neurons in 6 days (Scholz et al., 2011). We here describe functional properties of these cells as a fundament for the development of LUHMES-based pharmacological assays.

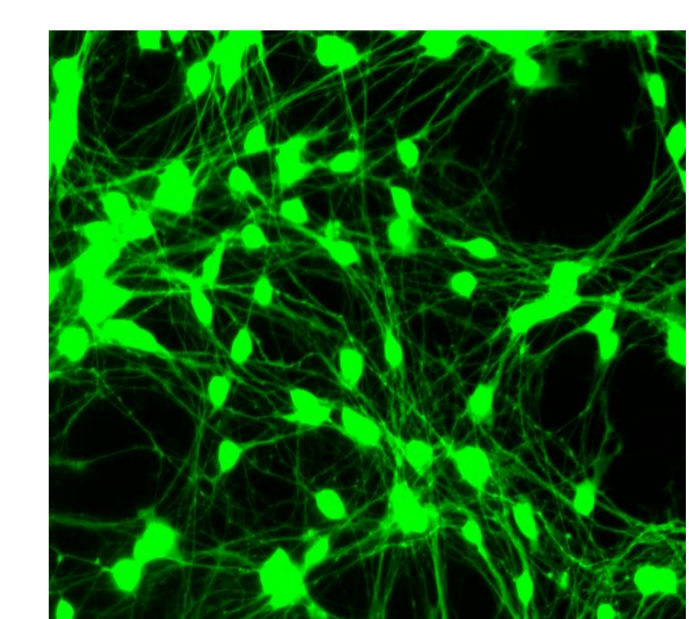
Materials & Methods

The LUHMES cells were cultivated for 9 days in differentiation medium in T75 or T175 flasks, coated with PLO (1 mg/ml) and Fibronectin (1 mg/ml). The recordings were performed with the patch clamp automat QPatch (Sophion Bioscience, DK) with 16X QPlates which enables parallel recordings of 16 independent experiments. For the experiments the cells were detached with 0.05% trypsin and resuspended in a concentration of 3-4 x 10⁶ cells per ml.



Materials & Methods

The LUHMES cells were differentiated for 2 days in T75 flasks, coated with PLO (1 mg/ml) and Fibronectin (1 mg/ml). The cells were further differentiated in a density of 20k per well in 384-well plates (Greiner Bio-One, GER), coated with 0.1% PEI. The Ca²⁺ imaging recordings were performed on day 9 of the differentiation with the HTS Functional Drug Screening System FDSS/μCELL (Hamamatsu Photonics, JP) combined with the Ca²⁺-sensitive dye Cal-520TM AM (AAT Bioquest, US) at 37°C.



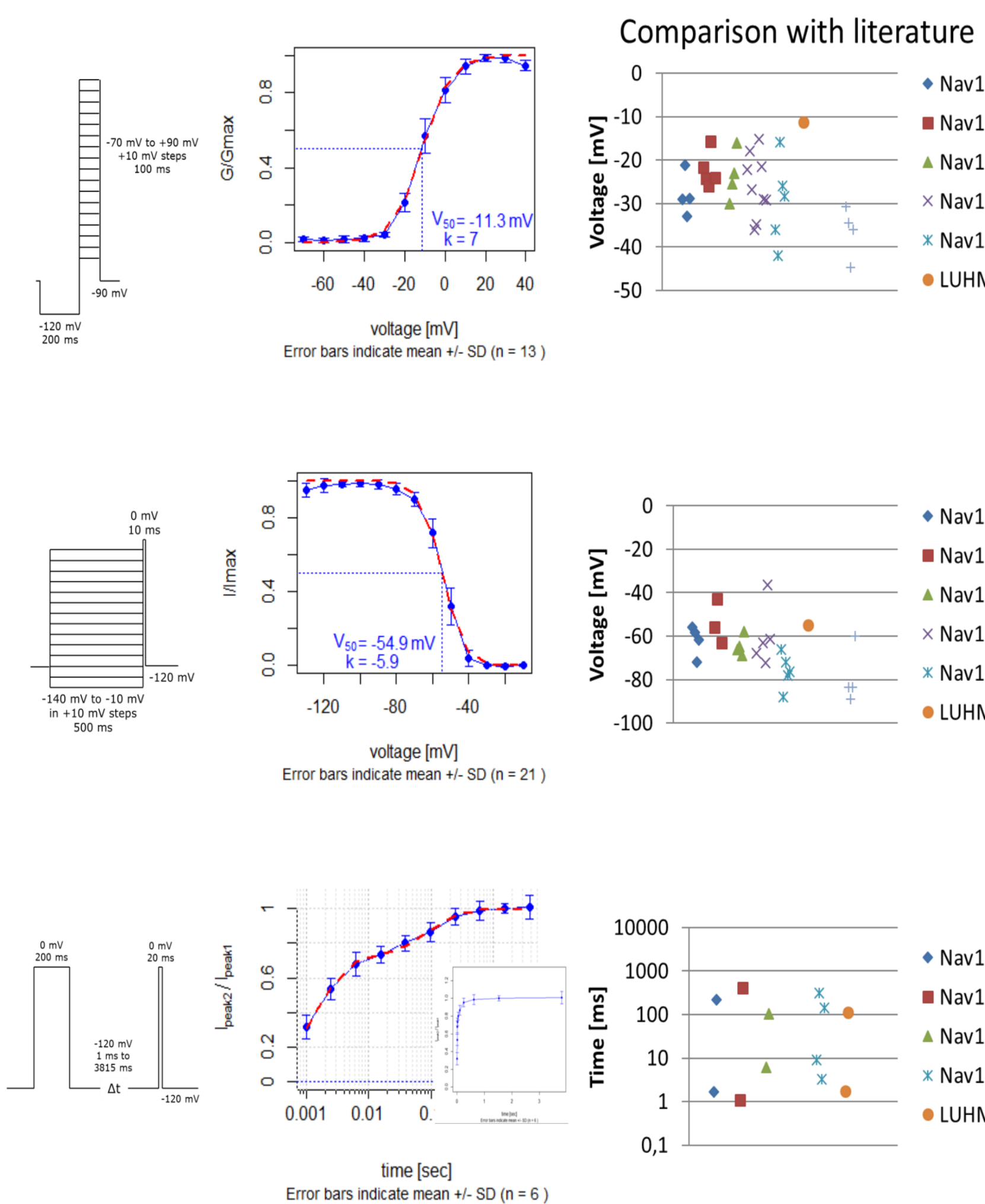
LUHMES cells stained with Cal-520TM on day 9 of the differentiation.



Results

Biophysical characterization of voltage-gated sodium (Na_v) channels using automated patch clamp

- Activation properties**
Evaluation of the normalized conductance (G/G_{max}) by estimating the half-maximal activation voltage (V₅₀) of -11.3 mV using a Boltzmann fit.
- Inactivation properties**
Analysis of the steady-state inactivation of the Na_v channels by calculating I/I_{max}. The Boltzmann fit determines a half-inactivation voltage (V₅₀) of -54.9 mV.
- Recovery from inactivation**
A bi-exponential fit was used to describe the obtained data. This resulted in two time constants of 1.71 ms & 112 ms.

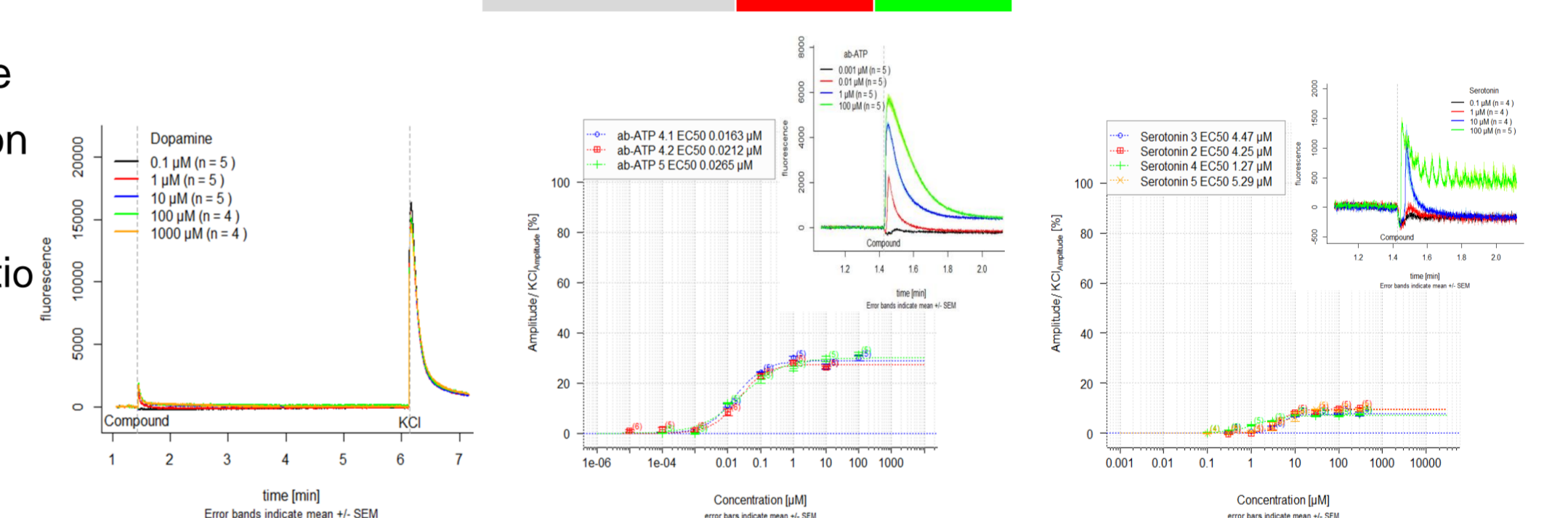


Results

Characterization of neurotransmitter receptors using HTS Ca²⁺ imaging

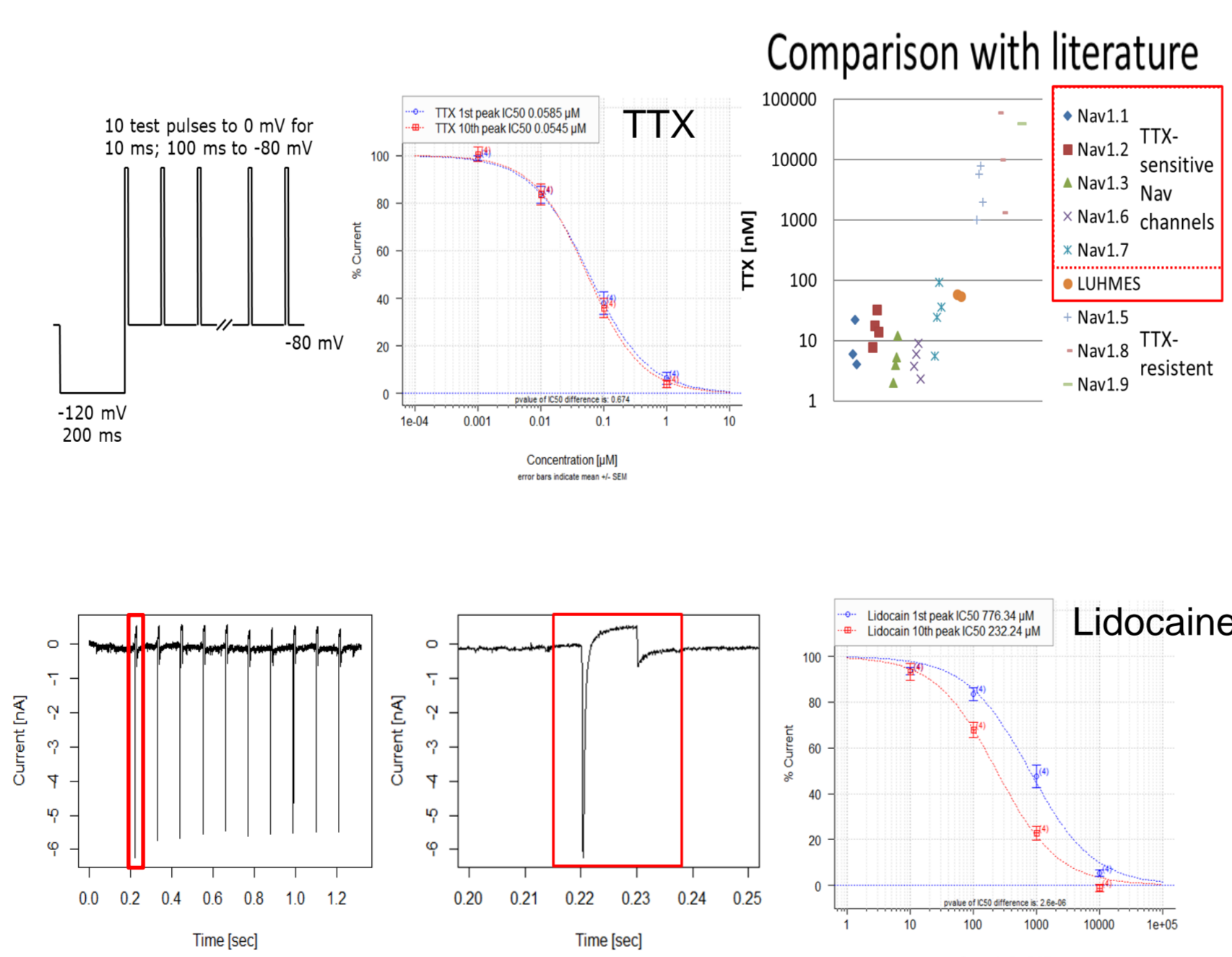
- Concentration-dependent effects of agonists of different neurotransmitter receptor were examined on various differentiations. The EC₅₀ values show a high consistency over several differentiations.
- Recording protocol:
 - 1.5 min control phase
 - Compound application
 - 4.5 min incubation
 - KCl (30 mM) applicatio

Receptor	Ca ²⁺ imaging analysis	mRNA
Dopamine	Green	Green
Serotonin	Green	Green
GABA	Green	Green
Glutamate	Green	Green
Nicotinic ACh	Green	Green
Muscarinic ACh	Red	Red
Adrenergic	Green	Green
Glycine	Green	Green
Histamine	Red	Red
Purinergeric	Green	Green
Adenosine	Red	Red



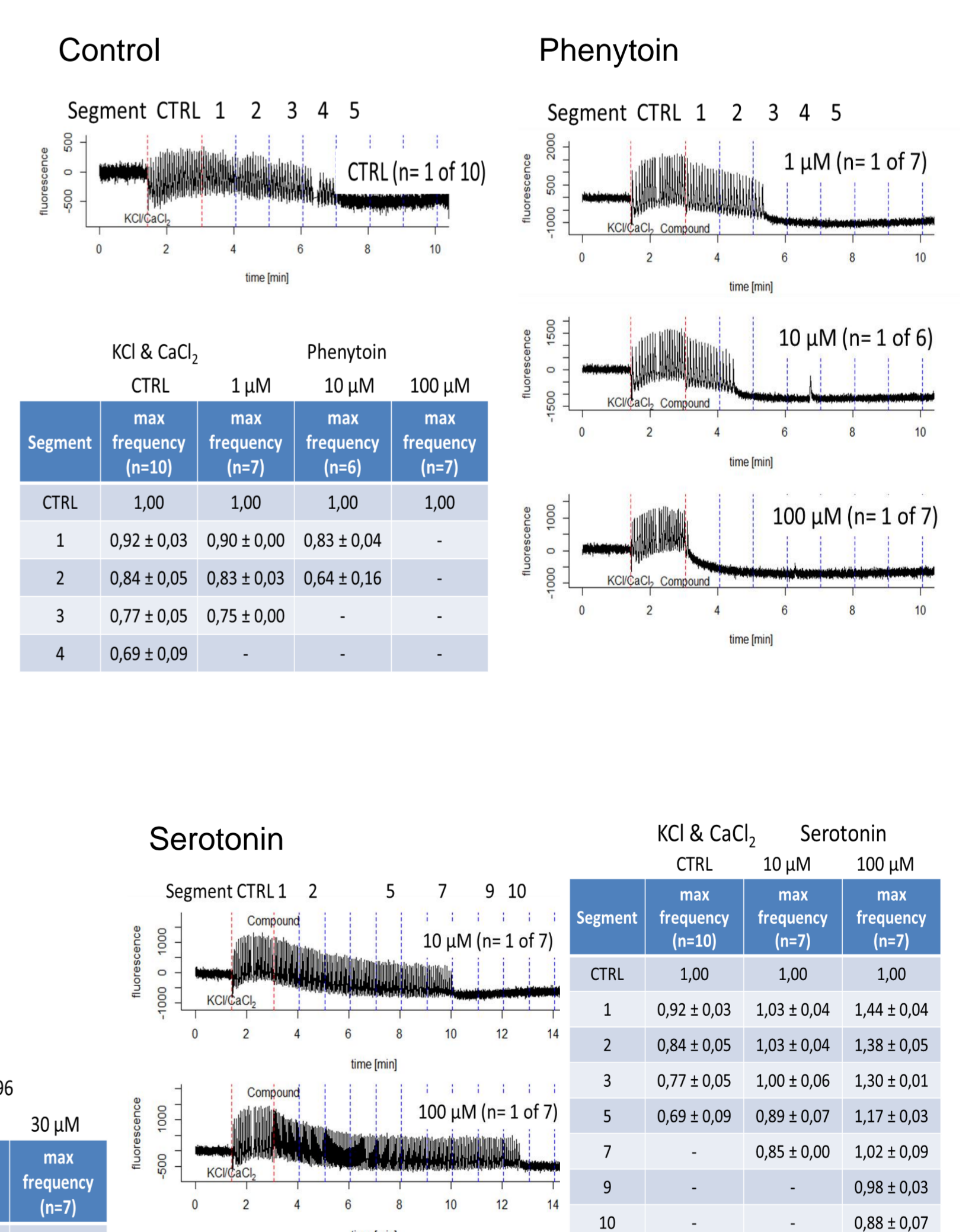
Pharmacological characterization of Na_v channels using automated patch clamp

- Effect of Tetrodotoxin (TTX) on Na_v channel currents to distinguish TTX-sensitive and TTX-resistant Na_v channels.** The IC₅₀ values suggest that only TTX-sensitive Na_v channel are present and TTX shows no use-dependent effect.
- Effect of Lidocaine (local anesthetic) which is a known use-dependent Na_v channel inhibitor.** The IC₅₀ values of the first and the last peak are significantly different.



Investigated of neuronal network activity using HTS Ca²⁺ imaging

- The oscillation frequency of the Ca²⁺ oscillations was examined with a FFT-based analysis.
- Hyperpolarization induced Ca²⁺ oscillations due to a synchronous activity of the neuronal network. Phenytoin (anticonvulsant drug) blocks Na_v channel currents and exhibits a concentration-dependent inhibition of the Ca²⁺ oscillations. Serotonin increased the oscillation duration and frequency.
- Serotonin (100 μM) induced Ca²⁺ oscillations were blocked by NNC 55-0396 (T-type Ca²⁺ channel blocker).



Conclusion

The results show that we were able to differentiate the cells derived from LUHMES cells with neuronal electrophysiological characteristics with low batch-to-batch variations. Addressing these neurons with the calcium imaging system could offer a great opportunity for a high-throughput assessment of the neurotoxic potential of novel drug candidates on a neuronal network in the future.

Acknowledgment for funding:

