

Preclinical Cytotoxicity Investigations in Stem Cells for Primary and Secondary Assays with an “All Inclusive” Approach

Corina T. Bot¹, Sonja Stölzle-Feix², Ulrich Thomas², Krisztina Juhasz², Leo Doerr², Matthias Beckler², Claudia Haarmann², Alison Obergrussberger², Markus Rapedius², Tom Götze², Marius Vogel², Susanne M. Rehnelt³, Tobias Brueggemann³, Philipp Sasse³, Michael George², Andrea Brüggemann², Rodolfo Haedo¹, Niels Fertig²

¹Nanon Technologies Inc., 1 Naylor Place, Livingston, NJ, 07039, USA

²Nanon Technologies, Gabrielenstr. 9, 80636 Munich, Germany

³Inst. für Physiologie I, Univ. Bonn, Life and Brain Center, Sigmund-Freud-Str. 25, 53127 Bonn

Human stem cell-derived cardiomyocytes (hiPSC-CMs) have recently proven to recapitulate key features of human cardiac electrophysiology in vitro.

Chip-based approaches allow parallel patch clamp recordings without compromising data quality or technical sophistication. We present high-throughput voltage and current clamp recordings of CiPA reference compounds. Since drug efficacies may vary with temperature, we present recordings at room and at physiological temperatures.

In addition to patch-clamping experiments, we present hybrid impedance (cell contractility) with MEA-like extracellular field potential (EFP) recordings. Experiments were complemented with optical stimulation of monolayers of hiPSC-CMs expressing the light-gated cation channel Channelrhodopsin2. This approach allows frequency-dependent drug screening and detection of potential side effects on Na⁺-, Ca²⁺- and repolarizing K⁺ channels. Furthermore, investigations of potential breaks in excitation-contraction coupling can accompany the ion channel screening with the final aim to enable a reliable automatized cardiac toxicity screening platform. As a proof-of-principle, pharmacological effects of CiPA compounds from each risk category will be presented.

Cytotoxic responses of cell monolayers involve metabolic or biochemical changes that affect the morphology of the cells, or reduce their overall viability. In that regard, effects of a number of reference compounds tested for long-term cytotoxicity in hepatocyte-like cells (i.e. Paracetamol), will be presented.