

# Complementary automated patch clamp, extracellular field potential and impedance recordings of iPSCs: safety screening tool box for the future

Elena Dragicevic<sup>1</sup>, Nadine Becker<sup>1</sup>, Sonja Stölzle-Feix<sup>1</sup>, Ulrich Thomas<sup>1</sup>, Krisztina Juhasz<sup>1</sup>, Leo Doerr<sup>1</sup>, Tom Goetze<sup>1</sup>, Markus Rapedius<sup>1</sup>, Daniel Millard<sup>2</sup>, Stacie Chvatal<sup>2</sup>, Anthony Nicolini<sup>2</sup>, Matthias Becker<sup>1</sup>, Michael George<sup>1</sup>, Andrea Brüggemann<sup>1</sup>, Niels Fertig<sup>1</sup>

<sup>1</sup>Nanon Technologies GmbH, Gabrielenstr. 9, 80636 Munich, Germany  
<sup>2</sup>Axion BioSystems, Inc., 1819 Peachtree Road, Suite 350, Atlanta, GA, 30309



## Patchliner



## CardioExcyte 96



## Axion Maestro

