

HIGH-END INSTRUMENTATION PROPOSAL

SyncroPatch 384i – Advance Ion Channel Screening

nanji[on]

Illawarra Health and Medical Research Institute (IHMRI)

Project Summary

The Illawarra Health and Medical Research Institute (IHMRI) is a not-for-profit charity that supports health and medical researchers in the Illawarra-Shoalhaven region of New South Wales, Australia. The aim of the research institute is to bring scientists, doctors and health professionals together to solve the greatest health challenges through research. IHMRI researches into better ways to prevent and treat disease and illness such as cancer, diabetes, mental illness and dementia.

The **SyncroPatch 384PE** is housed within the IHMRI's Electrophysiology Facility for Cell Phenotyping and Drug Discovery (E-Phys core). The instrument is used to measure electrical activity in up to 384 single living cells at a time, and their responses to different drugs. This speeds up the process of cell analysis and significantly reduces laboratory times for researchers.



SyncroPatch 384PE.
**Dr Alan Finkel AO (Chief Scientists, Australia) and
Professor David J. Adams (IHMRI)**

PROJECT SUMMARY

Since November 2018 [IHMRI's Electrophysiology Facility for Cell Phenotyping and Drug Discovery](#) has been successfully functionally characterizing multiple cell types and screening compounds on native and recombinant ion channels using the **SyncroPatch 384PE** and the PatchLiner octo.

We have successfully recorded native and recombinant voltage- and ligand- responsive currents in various stem cell derived models, neuroblastoma/glioma cell lines, primary and immortalized immune cells, as well as many ion channels stably expressed in various cellular backgrounds.

The mechanical and electrical stability of the system allows for precise and fast application of compounds making Nanion's automated patch clamp instruments invaluable tools for us.

Importantly, the team at Nanion provides outstanding customer support that is fast, knowledgeable and friendly.



SyncroPatch 384PE.
This Promotion is sponsored by Nanion Technologies GmbH, Ganghoferstraße 70a, 80339 Munich, Germany.

Please send an e-mail
SyncroPatch_Grant@nanion.de
for more information.

Nanion Technologies GmbH
Ganghoferstraße 70a
80339 Munich
Germany
www.nanion.de