

HIGH-END INSTRUMENTATION PROPOSAL

nan[i]on

SyncroPatch 384i – Advance Ion Channel Screening

Application Instructions

Title: Nanion Technologies GmbH, Grant Application "SyncroPatch 384i Grant: Advancing Ion Channel Assays"

By submitting the Application, you agree on the following rules:

- All entries must be submitted until the 30th of June 2020
- The declared winning Applicant of the Nanion Research Grant agrees to report work produced, data or results from the research proposal described in the Nanion Research Grant Application.
- If the winning Applicant plans to publicly present or publish the data and results of the research proposal described in the Application, the winning Applicant is required to inform Nanion Technologies GmbH prior to such publication or presentation.
- Nanion Technologies GmbH reserves the right to utilize the data derived from the winning Application for exhibition, promotion and marketing purposes in any medium.

Grant Committee and selection process

The Grant committee comprises the SyncroPatch 384i Technology specialists Dr. Andrea Brüggemann and Dr. Alison Obergrussberger from Nanion Technologies, GmbH, Germany and Rodolfo Haedo from Nanion Technologies Inc., USA.

The Grant Committee will determine the winning recipient based on (i) the best in-depth, novel application demonstrated with the SyncroPatch 384i platform; (ii) scientific merit of the application.

The decision of the Grant Committee is subjective, at the Sponsor's sole discretion and will be final and cannot be appealed.

Award Notification and Award

The winning application will receive a SyncroPatch 384i platform, 125 x 384-well consumables, training during the installation and consultation up to 9 months.

The winner may be posted by Nanion Technologies in social media channels and on the webpage.

Sponsorship

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