

Nanion Technologies Opens New Site in Copenhagen

Munich, Germany, June 28, 2016; Nanion Technologies is pleased to announce the opening of a new site in Copenhagen, Denmark. The site contains a fully equipped laboratory and state-of-the-art ion channel high throughput screening platforms, including the SyncroPatch 384/768PE. The new facility within Copenhagen University opens its doors for scientific collaborations, assay development support and product demonstrations.

Nanion is one of the leading providers of high-throughput automated patch clamp systems for industrial screening and academic research. With the inauguration of the facilities at Copenhagen University, and the recent recruitment of Søren Friis, Nanion is optimally positioned to offer excellent customer support to the constantly increasing customer base of the ion channel HTS-platform, the SyncroPatch 384/768PE. In addition, at the new site, Nanion will actively engage in scientific collaborations with researchers in pharma and academia.

Global Sales Director Frank Henrichsen says:

“With the rapidly growing customer base, and thereby increased demand for assay and technical support, we’re very happy to have Søren Friis on the team. His solid knowledge of ion channel applications and assay development combined with his dedication to customer support makes him an excellent fit for Nanion. The new facilities at Copenhagen University, and researchers in the area, such as Dr. Kirstine Callø and Dr. Dan Klærke, Faculty of Health and Medical Sciences, UCPH (Frederiksberg Campus) hold great promise for unique scientific collaborations and exciting projects between Nanion and researchers in the pharmaceutical industry and academia.”

Søren Friis, Senior Scientist continues:

“I am happy to join the Nanion team, and for the unique opportunity to build a lab in the midst of a flourishing academic community! There has always been a mutual appeal between Nanion and the scientific community, primarily because of the broad range of experimental possibilities offered by Nanion’s products. With the new site in Copenhagen, scientists can apply to use the instruments for their specific projects, and for everyone interested in live demonstrations, the SyncroPatch 384/768PE is ready for action.”

The inauguration during the 10th FENS Forum of Neuroscience | Copenhagen, will take place on July 4th, 2016 at the premises at IKVH, 100 Dyrnlægevej, Frederiksberg, where Nanion will hold an opening ceremony.

About Nanion Technologies:

Nanion Technologies is a one-stop-shop for ion channel drug discovery and screening technologies as well as sophisticated research instrumentation. Nanion was founded in 2002 as a spin-off from the Center for Nanoscience (CeNS) of the University of Munich. Nanion’s team has developed and successfully established four generations of automated patch clamp instruments for sophisticated and high throughput applications in ion channel research and drug discovery (Port-a-Patch, Patchliner and SyncroPatch product families). Further product lines are for cardiotoxicity screening (CardioExcyte 96), for parallel bilayer recordings (Orbit 16), and for parallel membrane transporter protein recordings (SURFE2R). Since 2014, Nanion distributes Axion’s multi-electrode array (MEA) systems in Europe and China.

Contact details:

Dr. Niels Fertig, CEO, Phone: +49 89 2189 97972, Email: info@nanion.de

Frank Henrichsen, Director of Global Sales, Phone: +45 40 87 29 17, Email: frank.henrichsen@nanion.de

Web: www.nanion.de