

## Søren Friis appointed Director of Global Customer Relations at Nanion

**Munich, Germany, February 7, 2018; Nanion Technologies is pleased to announce that Søren Friis has been promoted to Director of Global Customer Relations at Nanion. Søren Friis was most recently Senior Application scientist at Nanion. He holds a very strong track record in developing advanced assays for automated patch clamp systems and for providing excellent support to customers world-wide in pharmaceutical companies and CROs.**

Since the introduction of the high-throughput ion channel screening platform, SyncroPatch 384/768PE in late 2013 the platform has been installed globally in numerous ion channel screening and research laboratories. With a rapidly increased installed base mainly in pharmaceutical companies and CROs Nanion has launched an initiative to intensify the focus on the specific needs and requirements of pharmaceutical companies.

As an important part of the intensified focus on pharmaceutical companies Nanion has decided to promote Søren Friis to a newly created role as Director of Global Customer Relations. Søren has a large network in the pharmaceutical industry from customer support positions, and he has been working intensively with product and assay development on several automated patch clamp platforms since these platforms emerged more than a decade ago. During his work he has obtained a unique knowledge that pharma companies will benefit from in his new role.



Frank Henrichsen, Director of Global Sales at Nanion Technologies says: "For years, Søren and I have worked together, and I am thrilled to have him on our team in the key position as Director of Global Customer Relations. As well as his great personality, Søren

brings a wealth of knowledge and creativity and a thoughtful understanding of customer support into his new role. That will undoubtedly prove highly valuable in helping our customers to achieve their goals and timelines."



Søren Friis continues: "By listening to our customers, understanding their needs, and providing them with effective solutions, we earn their loyalty. This is how we have built Nanion's outstanding reputation. Excellent support is a key differentiator in our business and I look forward to working

with my colleagues at Nanion to further develop a first-class support framework in a competitive market."

### About Nanion Technologies:

Nanion Technologies is a leading provider of instrumentation for ion channel drug discovery and screening. Founded in 2002, Nanion has grown over the last 16 years to a company with over 100 employees worldwide. With headquarters in Munich, Germany, Nanion has subsidiaries in the USA, Denmark, Japan and China, as well as distribution partners in seven other countries. The Nanion 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.

### Contact Details:

Frank Henrichsen, Director of Global Sales  
Phone: +49 89 2190 95-0, E-mail: [info@nanion.de](mailto:info@nanion.de)  
Web: [www.nanion.de](http://www.nanion.de)