

Nanion nominated for The German Industry's Innovation Award

Munich, Germany, Jan 20th, 2009; Nanion is again nominated for a prestigious innovation award, due its impressive product portfolio of automated patch clamp systems. More than 350 companies competed in this year's Innovation Award and Nanion is finalist in the category "Start-Up Companies".

The German Industry's Innovation Award, also the world's first innovation award, has since 1980 annually nominated and rewarded the nation's most important scientific and technical innovations. Nanion is one of the five remaining candidates for the award and the winner will be announced on January 24th, 2009.

"To remain a profitable and leading provider of advanced and high quality ion channel drug screening systems, we constantly seek new inventions and solutions that are attractive to industrial and academic institutions." says Michael George, CTO of Nanion.

Dr. Niels Fertig, CEO of Nanion, continues *"Our three product families for sophisticated cell analysis are being used in the development of new medicines as well as in academic research. With the latest introduction of our high throughput patch clamp system, the SyncroPatch 96, the screening of ion channel drugs is revolutionized in terms of time and cost efficiency. We are happy and honoured that our products and technology are receiving such recognition by the nomination for this award."*

The SyncroPatch 96 is Nanion's next generation screening platform, capable of highly parallel recordings from up to 96 cells at a time. It will be launched at the Biophysical Society's Annual Meeting in Boston, MA, USA, in the end of February, 2009. The SyncroPatch 96 is the first platform on the market to date that supports giga-seal recordings at a throughput of 5000 data points per day.

About Nanion:

Nanion Technologies GmbH is a German Private Limited Company and 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 globally established two highly successful automated patch clamp instruments as enabling tools for sophisticated and high throughput applications for ion channel research and drug discovery.

Nanion's instruments use planar patch clamp chips which replace the traditional glass pipette used in the technique of patch clamping. Nanion was nominated in 2007 for Germany's most prestigious innovation award the **Deutscher Zukunftspreis** (German Future Prize, Federal President's Award for Technology and Innovation).

Contact details:

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

About the German Industry's Innovation Award

Under the patronage of the federal minister of economics and technology, the most influential scientific and technical innovations are each year nominated for the prestigious German Industry's Innovation Award (Innovationspreis der deutschen Wirtschaft), the world's first innovation award. The 29th annual award ceremony will take place in Frankfurt a.M. on January 24th, 2009.

For more information about the award, please visit: www.innovationspreis.com