

Nanion Acquires the SURFE2R Technology for Transporter Protein Analysis

Munich, Germany, February 26, 2012; Nanion acquires the well-established SURFE2R technology, allowing state-of-the-art measurements of transporter protein activity and function. The SURFE2R technology is an excellent enhancement to Nanion's existing products which until now covered ion channels and nanopores.

Nanion Technologies is one of the leading providers of high quality automated patch clamp systems. Nanion's efficient and versatile products drastically improve the data output in patch clamp-based ion channel screening, safety testing and scientific research. Since 2003, Nanion has launched three automated patch clamp platforms, the Port-a-Patch (2003), the Patchliner (2006), the SyncroPatch 96 (2010), and recently introduced the Orbit 16 – a device for parallel bilayer recordings. Existing products target ion channels and nanopores in cell- and cell-free membranes for highly detailed information about kinetics, function, and compound pharmacology.

Dr. Niels Fertig, CEO of Nanion Technologies, continues:

"By acquiring the SURFE2R technology developed by IonGate / Scientific Devices Heidelberg, Nanion enters yet another market segment. It is an excellent addition to our product portfolio and a unique platform since it allows direct measurements of transporter activity and function. Further on, the platform has been validated with a wide variety of membrane preparations ranging from bacteria to native cells and cell lines. With this acquisition, Nanion becomes a one-stop-shop for functional membrane protein analysis, extending its product portfolio from screening instrumentation for ion channels and nanopores to now also include transporter proteins. "

The SURFE2R product line will be launched during the second half of 2012.

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 of Nanion, Phone: +49 89 2189 97972, Email: info@nanion.de, Web: www.nanion.de