

Nanion increases throughput and cuts costs with a new industrial 96-channel patch clamp screening robot

Munich, Germany, November 18th, 2008; Today, Nanion announces the late-stage development of a new automated patch clamp platform: the **SyncroPatch 96**. Developed to meet the throughput demands of industrial ion channel drug screening and safety profiling, and with a price-per-data-point compatible with screening standards, the **SyncroPatch 96** will offer the highest throughput in the market for high quality HTS-oriented ion channel screening.

Following the successful market introduction of two automated patch clamp devices, the Port-a-Patch (2004) and the Patchliner (2006), Nanion now introduces the **SyncroPatch 96**. Nanion's Patchliner and Port-a-Patch platforms enjoy great popularity in both academic and industrial settings and have received enthusiastic user feedback in customer surveys such as the HTStec report. Building on their success, the new **SyncroPatch 96** vastly increases throughput while reducing the cost per data point to a level compatible with industrial ion channel screening requirements.

"There is a gap between the demands in ion channel drug screening and the capability of the high quality automated patch clamp platforms currently available on the market. Pharmaceutical companies want higher throughput and lower cost per data point, whilst maintaining data quality. The SyncroPatch96 will fill this gap, by providing high throughput, high quality patch clamp recordings, at a low enough cost to keep screeners happy." says Dr. Niels Fertig, CEO of Nanion.

The **SyncroPatch 96** acquires simultaneous recordings from 96 individual cells in a well-plate format and allows for screening of both ligand- and voltage-gated ion channels. The platform supports giga-seal recordings, continuous recording during compound application and addition of multiple compounds to each of the 96 cells. The **SyncroPatch 96** will be launched in 2009.

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