

## Nanion opens China office in Beijing

**Munich, Germany, October 24<sup>th</sup>, 2011; Today, Nanion announces the opening of yet another branch. To meet the needs of the rapidly expanding number of customers in China, the dedicated sales- and support team in Beijing now opens the doors to Nanion Technologies China.**

Nanion successfully entered the Chinese market over five years ago, and the installed base of Patchliner and Port-a-Patch instruments has grown constantly ever since. To cover the increasing interest and activities within the fields of ion channel screening, scientific and contract research, the subsidiary Nanion Technologies China is now officially open.

Andy Di, General Manager, Nanion Technologies China, says:

*"Nanion's customer base in China has grown rapidly over the past couple of years. Nanion's products are very appreciated here because of their vast experimental freedom paired with increased throughput and affordable consumables. Because of their high quality and reliability, products like the SyncroPatch 96, the Patchliner and the Port-a-Patch are attractive for both industrial screening purposes and for dedicated academic research."*

Niels Fertig, CEO of Nanion Technologies GmbH, continues:

*"Because of the expanding Chinese market and customer base, it was a natural step to expand in China. With the new subsidiary we are able to cover the ever-increasing demand for product demonstrations and installations as well as to continue giving our Chinese customers premium support. We definitely feel that we have the right people in place for this important step in Nanion's history, and we look forward to a fruitful collaboration with the Nanion team in China. "*

Nanion Technologies China offers live product demonstrations and customer training in Beijing and Shanghai, or at the customer site.

### About Nanion

Nanion Technologies GmbH is a German, privately held, limited company, founded in 2002 as a spin-off from the Center for Nanoscience (CeNS) of the University of Munich, Germany. Nanion's team has developed and globally established three successful automated patch clamp instruments as enabling tools for sophisticated and high throughput applications for ion channel research and drug discovery.

Nanion's high quality instruments employ planar patch clamp chips which replace the traditional glass pipette used in the conventional patch clamp technique. In 2009 Nanion was awarded the German Founders Award (Gründerpreis), and the Innovation Prize Step Award. Next year, Nanion celebrates its 10<sup>th</sup> anniversary. Nanion has two branches: Nanion Technologies Inc., North Brunswick, NJ, USA, opened in 2009, and Nanion Technologies China, opened 2011.

### Contact details, Nanion Technologies GmbH:

Niels Fertig, CEO, Phone: +49 89 2189 97972, Email: [info@nanion.de](mailto:info@nanion.de), Web: [www.nanion.de](http://www.nanion.de)

### Contact details Nanion Technologies China:

Andy Di, General Manager, Phone: +86 1082176388 Email: [Andy.Di@nanion.de](mailto:Andy.Di@nanion.de) Web: [www.nanion.de](http://www.nanion.de)