

Nanion Technologies GmbH

Pettenkoferstr. 12
80336 Munich
Germany

The logo for Nanion Technologies GmbH, featuring the word "nanjion" in a white, lowercase, sans-serif font. The letters "j" and "i" are stylized with a vertical bar through them. The logo is set against a solid black rectangular background.

fon: 089 2180 75 260
fax: 089 2180 75 001

web: www.nanion.de
e-mail: info@nanion.de

contact: Dr. Niels Fertig (CEO)

Nanion receives Bavarian Innovation Award 2005

Munich, Germany – Octobre 2005 – The German nanobiotech company Nanion Technologies GmbH was recently awarded the Bavarian Innovation Award 2005 and is Bavarias company of the year (small/medium enterprise (SME)).

Nanion received the award for its chip-based patch clamp technique, which allows to perform sophisticated electrophysiological measurements from cells in an automated and parallel manner. Nanions entry level device, the Port-a-Patch[®], supersedes the common pipette based patch clamp technique in ease of use and throughput and doesn't require a microscope, vibration isolation or a micromanipulator. The Port-a-Patch[®] uses an all electrical read-out for automatically positioning and electrically contacting cells on the chip. With the Port-a-Patch[®], Nanion offers the world's smallest patch-clamp workstation. This innovative drug discovery technology makes patch clamp available also to non-electrophysiologists.

This innovation and its successful market introduction has been award the prestigious prize, which clearly is an important step and great commendation for Nanion.

Nanion

Nanion Technologies GmbH is a spin-off from the Center for Nanoscience (CeNS) of the University of Munich (LMU). Nanion combines bio- and nanotechnology in a company serving the life sciences industry by offering instrumentation for increasing the speed and efficiency of the drug discovery process in ion channel drug discovery. Nanion develops, produces and markets instrumentation and chip consumables for automated patch clamping. For more information, please contact us at info@nanion.de or visit our website at www.nanion.de.

The Port-a-Patch[®]

The Port-a-Patch[®] is a complete patch clamp setup with minimum foot print and low maintenance requirements. The system uses Nanion's planar patch clamp chips and provides high quality data. Great ease-of-use is accomplished by automated cell positioning and sealing. The Port-a-Patch[®] enables fast fluid exchange on the chip, suitable not only for voltage gated, but also for ligand gated ion channels. Even the exchange of intracellular solution is possible, enabling perfusion on both sides of the cell membrane. The Port-a-Patch[®] offered by Nanion is a valuable tool for target validation and safety pharmacology (hERG screening).