

Nanion Introduces Platform-Certified Cell Lines

Munich, Germany, November 08, 2012; Nanion's automated patch clamp platforms are known for their high cell compatibility, even for primary cells and stem cells. Nanion now introduces "Nanion Approved Cells", where commercially available cell lines from several providers have been evaluated and certified based on cell and platform performance. The approved cells can be directly purchased through Nanion, so that Nanion's customers get a complete package of APC platform and validated assay.

A large number of cell lines from several providers, including Millipore and CreaCell have been validated on Nanion's three automated patch clamp platforms, the Port-a-Patch, Patchliner, and the SyncroPatch 96. Parameters such as average seal-resistance, access resistance, cell recording longevity, ion channel expression and success rates for completed recordings were investigated and documented in a standardized way. Success rates for completed recordings were typically above 70%. With Nanion Approved Cells, the user can rest assured that the assay will run optimally, with minimal or no assay development required.

Atsushi Ohtsuki, Customer Support Manager, Nanion Technologies, Japan says:

"The Nanion Approved Cells will be much appreciated by Japanese customers, since the optimization work already has been done by Nanion Technologies, which saves time and reduces the risk of an investment. Additionally, it is also a guarantee, that the customers will encounter minimal problems when starting a new assay on Nanion's automated patch clamp platforms."

Dr. Niels Fertig, CEO of Nanion, continues:

"The certification of commercially available cell lines makes it easier for existing and new customers to invest in new cell lines and our platforms without having to worry about lengthy assay development. It is a well-known fact that Nanion's platforms show great compatibility with a broad range of cells and ion channels, which also is confirmed by the long list of Nanion Approved Cells and scientific publications in high impact journals such as Nature or PNAS. Nanion also certifies cell lines on-demand, meaning that if a customer has a particular cell line of interest, we evaluate and document the cell performance on our platforms, free of charge."

About Nanion Approved Cells

Nanion Approved Cells are commercially available cells, certified on Nanion's three automated patch clamp platforms. Cell and platforms performance as well as ion channel current expression is thoroughly investigated and summarized in a data sheet. The approved cells can be directly purchased from Nanion, so that Nanion's customers get a package of APC platform and validated assay. Nanion also does cell line certification on-demand, if the cell line of interest cannot be found in the list of approved cells. The full list of tested cell lines can be found on Nanion's website.

About Nanion Technologies:

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 three highly successful automated patch clamp instruments as enabling tools for sophisticated and high throughput applications for ion channel research and drug discovery.

Nanon'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).