

The SyncroPatch 384 is installed at the FDA for cardiac safety pharmacology studies

New Jersey, USA, March 10th 2022: Nanion is pleased to announce that a SyncroPatch 384 instrument has been installed at the Food and Drug Administration (FDA) at their Silver Spring site in Maryland. The SyncroPatch 384 will be used for high throughput automated patch clamp experiments within cardiac safety.

About the SyncroPatch 384

The SyncroPatch 384 is a high throughput automated patch clamp instrument for recording from up to 384 wells simultaneously.

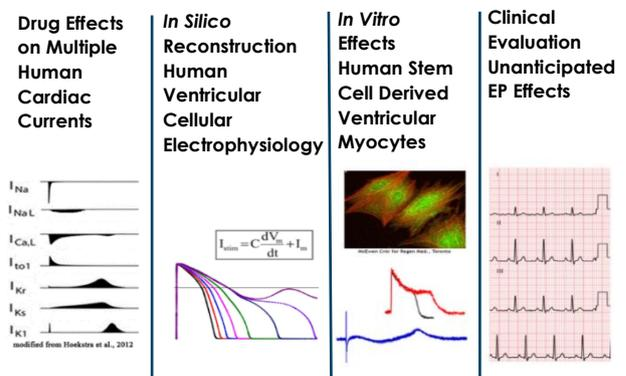


The system comes equipped with temperature control for heated experiments at physiological temperature, or cooled experiments below room temperature. It is highly flexible and easy-to-use and experiments can be performed using different solutions, with or without fluoride and high divalent solutions. The 32-well mode can be used for smaller screening projects and assay development.

About cardiac safety testing

The ICH S7B non-clinical guidance was introduced in November 2005 and required all new drugs to be tested for activity on I_{Kr} carried by hERG-expressed in recombinant cell lines using the patch-clamp technique, as blockade of the hERG channel has been attributed to drug-induced long QT syndrome and potentially fatal Torsade de Pointes (TdP). Since the introduction of these guidelines, very few drugs have been withdrawn from the market due to pro-arrhythmic complications. In 2013 the Comprehensive *in vitro* Proarrhythmia Assay (CiPA) initiative was started to further improve risk prediction, by accounting for drug block of cardiac ion channels (e.g., Nav1.5 and Cav1.2) that may mitigate the risk of hERG block. This effort is to ensure that drugs that reach the market are safe, by identifying candidate compounds with the potential to cause potentially fatal arrhythmia, but also to reduce the number of false positives ensuring that potentially useful (and safe) drugs do reach the market.

Comprehensive *in Vitro* Proarrhythmia Assay: Four Components



The four components of the CiPA initiative.



Chief Operating Officer at Nanion Technologies Inc., Rodolfo Haedo, was thrilled to announce the installation at the FDA saying, 'Cardiac safety testing of compounds is paramount to

ensuring safe compounds reach the market, and typically ion channel data submitted to the FDA are generated by manual patch clamp systems. The SyncroPatch 384 has now been installed at the FDA's Silver Spring site, and it will be used to evaluate how automated patch clamp data may be used in the regulatory context. Data generated by the SyncroPatch 384 will be compared with those concurrently generated using the manual patch clamp method, and we look forward to this comparison'.

About Nanion Technologies

Nanion Technologies is a leading provider of instrumentation for ion channel drug discovery and screening. Founded in 2002, Nanion has grown over the last 20 years to a company with over 100 employees worldwide. With headquarters in Munich, Germany, Nanion has subsidiaries in the USA, Japan, China and Denmark, as well as distribution partners in seven other countries. Nanion's team has developed and successfully established four generations of automated patch clamp instruments for sophisticated and high throughput applications in ion channel research and drug discovery (Port-a-Patch, Patchliner and SyncroPatch product families). Further product lines are available for cell monitoring and cardiotoxicity screening (CardioExcyte 96/FLEXcyte 96), parallel bilayer recordings (Orbit family), and parallel membrane transporter protein recordings (SURFE²R).

For more information, please visit:

www.nanion.de

Contact:

Rodolfo.haedo@naniontech.com



Nanion Technologies GmbH Director of Scientific Affairs and CiPA panel expert, Dr. Sonja Stölzle-Feix said, 'The CiPA initiative is an ongoing project aimed at ensuring cardiac safety of

new compounds, whilst at the same time making sure that safe drugs don't fail early on in the pipeline due to hERG block when they are not pro-arrhythmic. Both the FDA and Nanion Technologies, amongst other institutions and companies, have been heavily involved in standardizing experimental protocols for patch clamp experiments. We are looking forward to working with the FDA to ensure efficient transfer of their manual patch clamp assays onto the SyncroPatch 384.'

About the FDA

The Food and Drug Administration is responsible for protecting the public health by ensuring the safety, efficacy, and security of human and veterinary drugs, biological products, and medical devices; and by ensuring the safety of our nation's food supply, cosmetics, and products that emit radiation.

FDA also has responsibility for regulating the manufacturing, marketing, and distribution of tobacco products to protect the public health and to reduce tobacco use by minors.

FDA is responsible for advancing the public health by helping to speed innovations that make medical products more effective, safer, and more affordable and by helping the public get the accurate, science-based information they need to use medical products and foods to maintain and improve their health.

FDA also plays a significant role in the Nation's counterterrorism capability. FDA fulfills this responsibility by ensuring the security of the food supply and by fostering development of medical products to respond to deliberate and naturally emerging public health threats.