

Nanion Technologies Opens New Office in Silicon Valley

Newark, USA, August 24, 2018; Nanion extends its reach by opening a new US office in the San Francisco Bay Area, California. In order to support the increasing demand of Nanion's products and to continue providing premium support and customer service, Nanion now officially expands its physical presence on the West Coast.

Nanion opens the new office in a state-of-the-art incubator site strategically located in Silicon Valley where several innovative start-ups and small companies work together in a vibrant ecosystem of collaboration. Dr. James Costantin, Director of Customer Relations, opened the office in July of 2018 and we are welcoming new scientific collaborations, as well as providing assay support for both our pharmaceutical and academic customers. A newly hired Field Application Scientist will also be located out of the Silicon Valley office.



Rodolfo Haedo Senior Vice President North America, says:



"I am delighted to see Nanion continue to grow so rapidly in the US. As we expand to the West Coast, we can further strengthen the focused and attentive customer support that has helped build our success. As the sales of our SyncroPatch 384/768PE HTS systems have substantially increased over the last years, we have decided to expand both our infrastructure and personnel in this region full of young energetic startup companies. In addition to the office and new West Coast Field Application Scientist position, we anticipate future employment opportunities in order to help support and foster the rapidly increasing customer base. We are all very much looking forward to this new phase we are entering at Nanion."

Dr. James Costantin, Director of Customer Relations, continues:



"The proximity of Nanion's West Coast office to the rapidly expanding presence of pharmaceutical companies heavily investing in both the South San Francisco and San Diego regions makes the West Coast office ideally situated to serve our customers. In addition to providing office space to serve our customers, the new location allows Nanion to provide rapidly responding Application Support across all US time zones. We are driven to ensure that our customers succeed at all levels of their ion channel Drug Discovery endeavors, including Assay Development, Screening Campaigns, SAR/Medicinal Chemistry, and Counter Screening. Our customer's success is extremely important to us, as it is how we measure our success."

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 16 years to a company with over 100 employees worldwide. With headquarters in Munich, Germany, Nanion has subsidiaries in the USA, Denmark, Japan and China, 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 for cardiotoxicity screening (CardioExcyte 96), for parallel bilayer recordings (Orbit family), and for parallel membrane transporter protein recordings (SURFE²R).

Contact Details:

Rodolfo Haedo, SVP, Phone: (862) 221-2790
Email: rodolfo.haedo@naniontech.com

Dr. James Costantin, Director of Customer Relations,
Phone: (862) 596-9408

Email: james.costantin@naniontech.com

Web: www.naniontech.com

39655 Eureka Drive, Newark, CA 94560-4806, USA