

An aerial photograph of a person surfing a massive, curling wave. The water is a deep, dark blue, and the wave's crest is white with foam. The surfer is a small figure in the lower center, riding the base of the wave. The overall scene is dynamic and powerful.

SyncroPatch 384i

When performance matters.

#automatedpatchclamp

nanji[on

reliability scalability flexibility

*“The field of ion channel screening is challenged by an increasing demand for high throughput by Big Pharma and CROs, who need to **reliably screen** more compounds on diverse targets.*

*On the other hand, researchers have smaller, more diverse development projects that need a **higher flexibility**. The SyncroPatch 384i is ideally suited for both purposes.”*

*Dr. Niels Fertig, PhD, Co-Founder & CEO
Nanion Technologies*

Your new SyncroPatch 384i

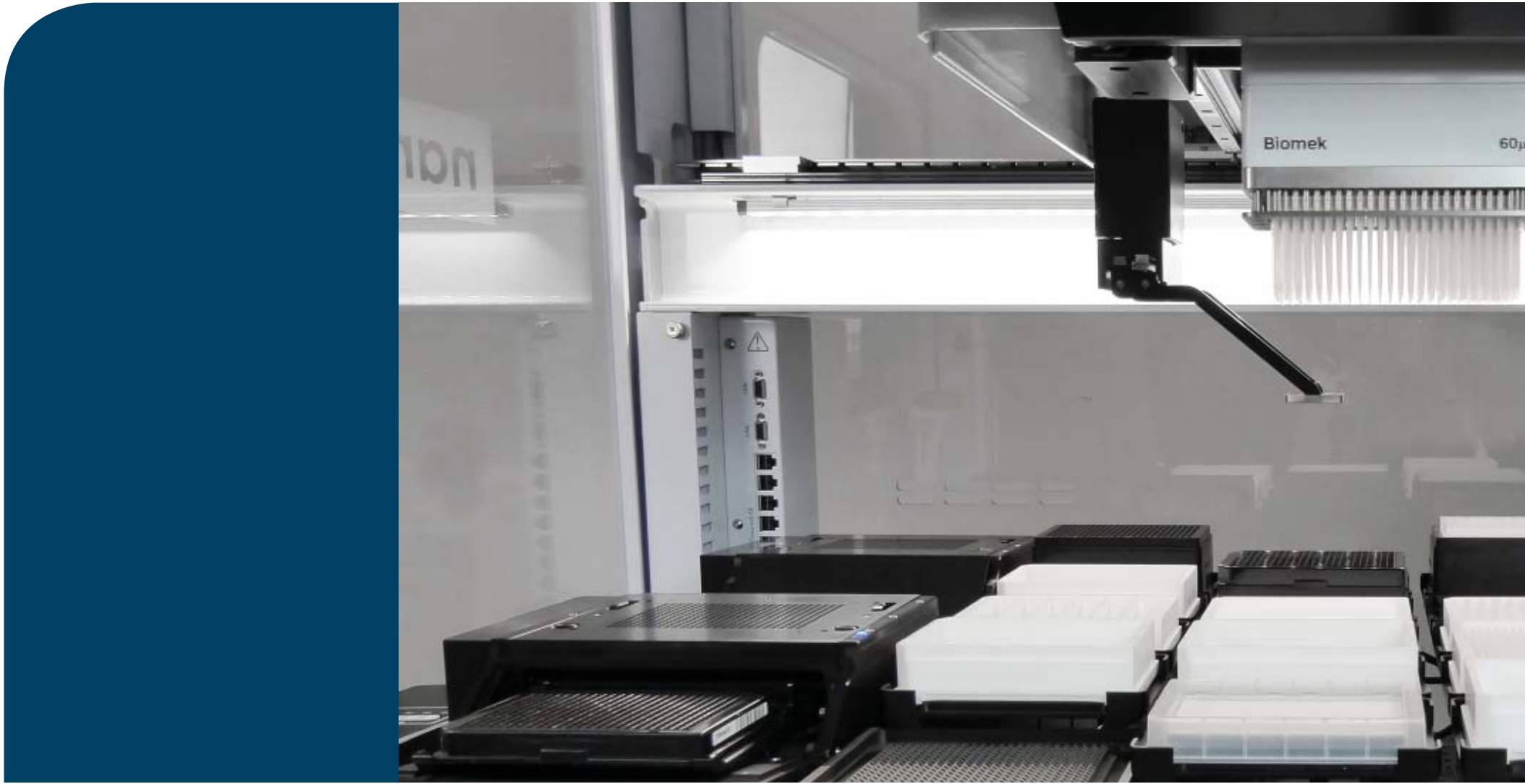
The best of both worlds.

Starting a thrilling new era

Fifteen years ago we introduced our first automated patch clamp system. During that time, our portfolio and our technology has constantly improved, as we have been listening to the needs of our customers in order to offer the best possible innovative instruments for APC and beyond.

The new SyncroPatch 384i provides effortless screening coupled with high flexibility and reliability, due to the combination of our patch clamp module with Beckman Coulter's new i-Series liquid handling robot.





Powered by Biomek i5

Designed for your workflow priorities.



Welcome to the new liquid handling

- Large deck space
- Linear motion control
- Fast pipette dispense speed
- Independent movement of gripper and pipetting head
- Selective tip pipetting
- LED-illuminated deck

*Enabling
innovation*



384 PS

"Pferdestärke" (PS or horsepower) is a "unit of measurement of power, or the rate at which work is done".

The SyncroPatch 384i contains a 384 "PS" module for 384 recordings in parallel. The new SyncroPatch 384i builds on the success of the SyncroPatch 384PE, which has been globally established as the preferred automated patch clamp workhorse in Pharma, Biotech, CRO and academia, whilst further improving flexibility, ease-of-use and reliability.

- 384 channel automated patch clamp system
- Suitable for use in all phases of drug discovery
- Optional add-ons: Temperature control; current clamp; second module for doubling your throughput to achieve up to 40.000 dp/day

Your Software Cockpit

Even more powerful than before

PatchControl and DataControl

Editing and analyzing the easy way

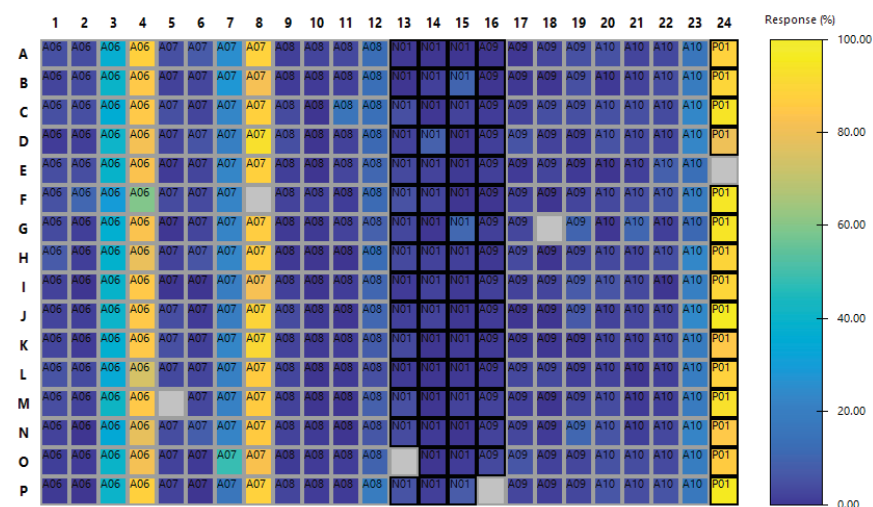
New tools

- Z-prime calculation
- Heat maps

Enhanced ease-of-use

- True partial plates
- Method Launcher
- Guided Labware Setup
- Windows 10 OS

Heat maps



Get a quick overview - e.g. response (%) to compounds across 384 wells.

Maximizing your Data Output

Your goals and everyday challenges

We maximize instrument uptime and data output at every stage

You'll never walk alone: Get supported by our 25 highly skilled electrophysiologists – with a total of more than 400 man-years of experience, you will receive unsurpassed application and technical support.

Benefit from our fast service and system diagnosis via onboard cameras.

Choose between two service plans: Premium and Basic. We are your single point of contact for the entire system.

Improved Simplicity You will love

You have intuitive software tools at hand for setting up your experiment with the Guided Labware Setup. The new Method Launcher allows for easier operation for non-experts. With the on-board cameras you gain control throughout your experiment.

- Progress Bar with Milestones
- On-deck compound plate preparation
- 360 ° multi-color status light bars for system status
- Integration into automated HTS environments
- High experimental flexibility - experiments can be modified on the fly
- Economical cost per data point
- Improved assay development - option for partial plate measurements (down to 1/12 of a measurement plate)
- Optimized handling of sticky compounds
- Internal perfusion during recording



Nina and Marius, just two of our experts who are happy to introduce you to our new SyncroPatch 384i.

**schedule your
demo now!**

<https://www.nanion.de/en/schedule-a-demo.html>



*“With the SyncroPatch 384i we are **improving the user experience** by offering a simpler, more efficient and more reliable HTS automated patch clamp system.”*

*Dr. Claudia Haarmann, Product Manager of SyncroPatch 384i and
Nanion’s Director of Product Implementation (APC)*

improving
the user
experience

Nanion Europe

info@nanion.de

phone: +49 89 2 190 95-0

www.nanion.de

Nanion USA

info@naniontech.com

phone: 1-888-9-NANION

www.naniontech.com

Nanion China

andy.di@nanion.cn

phone: +86 10 82 17 6386

www.nanion.cn

Nanion Japan

ohtsuki@nanion.jp

phone: +81 3 6457 8776

www.nanion.jp