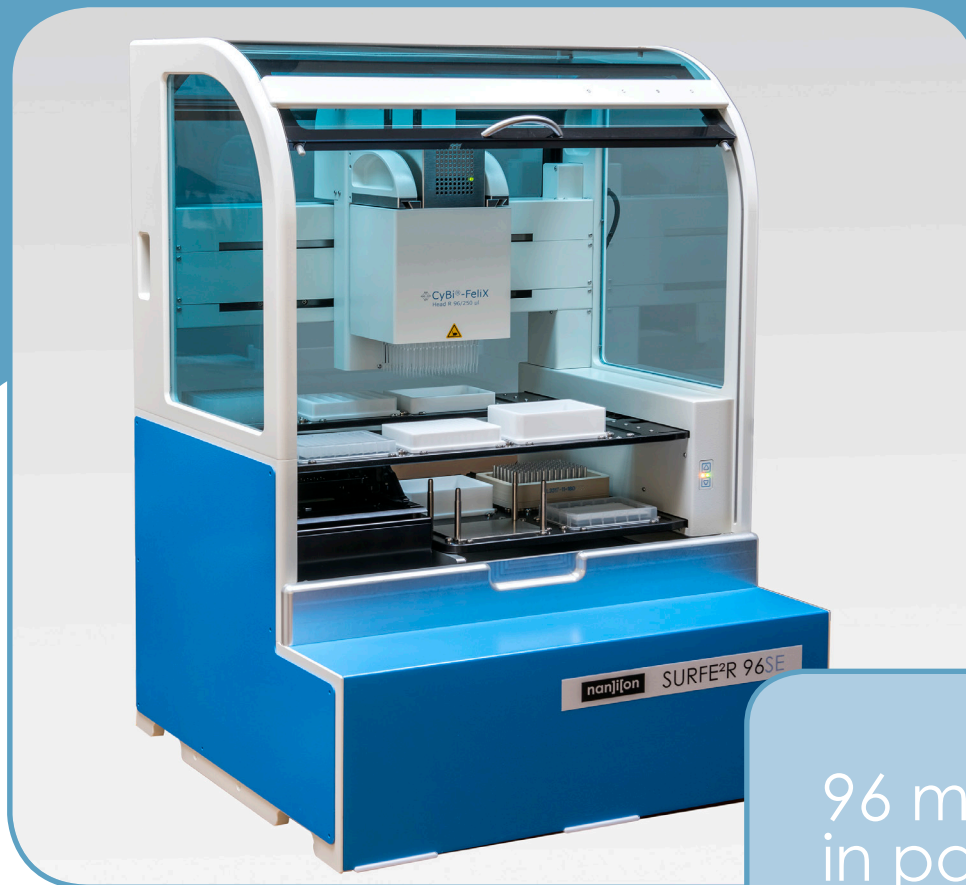


SURFE²R 96SE

Official launch of HTS device for electrogenic transporters and pumps



This is
New:

Visit us at booth #405
to see our new
SURFE²R 96SE

The new SURFE²R 96SE is the first high throughput device for the electrophysiological analysis of transporters and pumps

The new SURFE²R 96SE offers you:

- 96 datapoints in less than 5 minutes
- Label-free real-time measurements
- No radioactivity
- An ultrafast solution exchange with a time resolution down to 5 ms
- A superior signal to noise ratio

96 measurements
in parallel

Based on SSM
technology

Robust and
reliable

nanjion

Nanon Europe
info@nanion.de
Phone: +49 89 2 18 99 79 72
www.nanion.de

Nanon USA
info@naniontech.com
Phone: 1-888-9-NANION
www.naniontech.com

Nanon China
info@nanion.cn
Phone: +86 10 82 17 63 86
www.nanion.cn

nanjion

The new SURFE²R 96: HTS of transporter and pumps

"The SURFE²R 96SE is a full parallel 96 well-based system, generating 96 data points in less than 5 minutes. The device is designed for activity measurement of electrogenic membrane transporters and pumps. It focuses on targets not accessible for patch-clamp, such as low turnover targets and targets localized in intracellular membranes or membranes of small cells like bacteria. Solid Supported Membrane (SSM) based electrophysiology enables characterization of transported substrates, inhibitors, modulators and cofactors and is used to derive kinetic parameters such as EC_{50} , IC_{50} , rate constants and more. The measurement uses a rapid solution ex-

change with a time resolution down to 5 ms to even characterize fast binding kinetics using arbitrary substrates. The fact that membrane fragments or liposomes are measured instead of living cells yields in a high reproductivity and robustness of this technology and a superior signal to noise ratio of the electrical measurement. We are very excited to introduce this innovative high throughput system. With this, we can offer our clients a screening platform for these difficult targets in a robust manner and meaningful throughput"

Dr. Maria Barthmes, Product Manager of the SURFE²R product family, Nanion Technologies GmbH

Do not miss:

- To learn if the SURFE²R 96SE is the perfect solution for your transporter, pump or ion channel project.

Visit our booth # 405 and talk to our SURFE²R specialist:

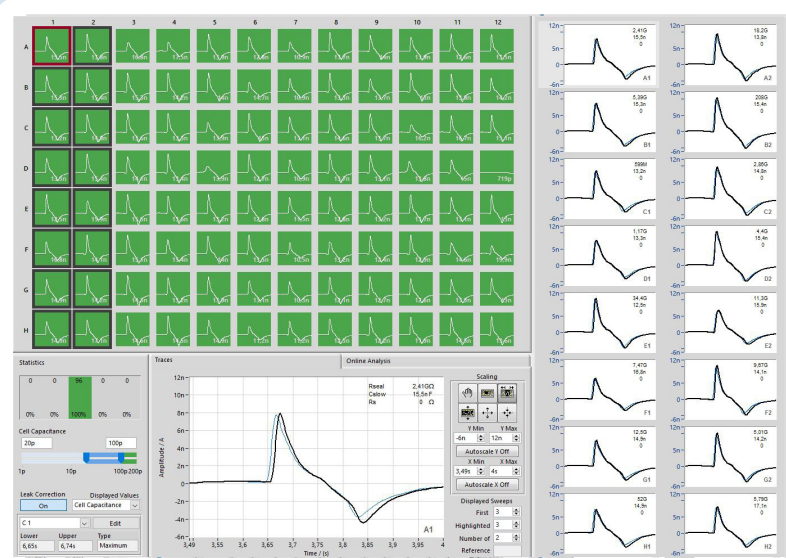
- Maria Barthmes, product manager of the SURFE²R product family

Visit our Exhibitor presentation (Tuesday, 12:30 PM, room 221):

- "Measure More Membrane: Cells and Bilayers on the Port-a-Patch, SURFE²R, Orbit 16 and Orbit Mini"

Visit our Poster presentation (Monday 2:45 PM, 1348-Pos, B416)

- "Functional characterization of the SLC-Transporter PEPT1 and OCT2 by electrophysiological real-time measurements using a high throughput system"



Activation of PepT1 on SURFE²R 96SE:

Graphical user interface of the screening software of the SURFE²R 96SE: Traces of PepT1 recorded from membrane preparations of PepT1 overexpressing CHO cells are presented. IC_{50} determination with five concentrations, duration 15 minutes. The PepT1 signal is evoked by 20 mM glycylglycine.