

Port-a-Patch

New functions and new design



This is
New:

New design

New Temperature
control

New External
perfusion

Visit us at booth #405
to see our new
Port-a-Patch

The new Port-a-Patch

- Proven reliability in new design
- Still smallest patch clamp rig in the world
- Heatable AND coolable temperature control

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



The New Port-a-Patch

"The Port-a-Patch is a miniaturized and easy-to-use patch clamp system that is well established and accepted among CRO, pharma and academia since its market entry 12 ago. Here we introduce the new Port-a-Patch reloaded with a new design and a completely newly developed external perfusion system. The new perfusion system has an integrated temperature control enabling you to perform experiments at higher AND lower temperatures in regard to room temperature. In addition more add-ons like internal perfusion, microscope slide and Low Noise Unit are available.

Using the Port-a-Patch is straightforward and easy –the user simply adds solutions and cells onto the disposable recording chip, where a cell is automatically captured and sealed by suction using a computer controlled pump.

The Port-a-Patch system is comprised of the Port-a-Patch recording unit, a Suction Control unit, a HEKA EP-C10USB amplifier and a computer "

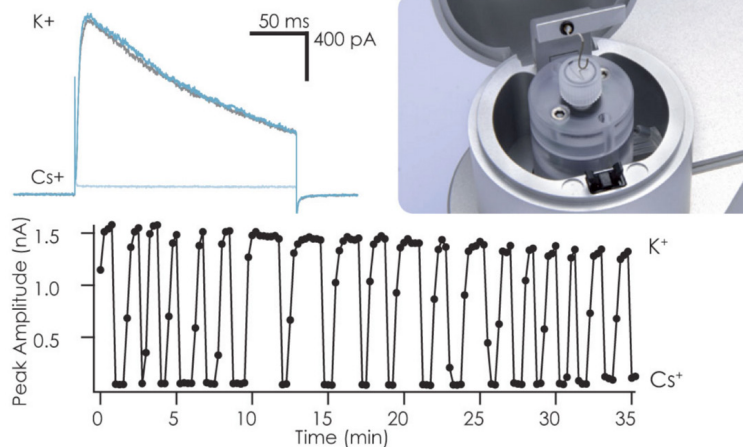
Dr. Patrick Mumm, Product Manager Port-a-Patch, Nanion Technologies GmbH

Do not miss:

- To learn about the new design of the Port-a-Patch
- To learn about the new features of the temperature control:
Cooling and heating is especially interesting for the measurement of temperature sensitive TRP channels.
- To learn about the internal and external perfusion capabilities of the Port-a-Patch
- To find out how simple it is to use the Port-a-Patch for your studies

Visit our booth # 405 and talk to our Port-a-Patch specialists:

- Andrea Brüggemann, George Okeyo, Corina Bot, Niels Fertig



Continuous Internal Perfusion

$K_v1.3$ currents (blue), endogenously expressed in Jurkat cells, were rapidly blocked by internal perfusion of Cs^+ (light blue), and fully recovered after washout with K^+ (grey). Internal solution replacement was repeated 19 times and the recording was stable for over 35 minutes, as shown in the lower graph.