

Suction Control Pro

– keeping up the pressure

- **Pressure clamp extension for Nanion's second generation Suction Control**
- **Improved stability and lifetime**
- **Fast and easy auto-calibration and self-test**
- **Faster pressure regulation**

The second generation of the Suction Control unit for the Port-a-Patch® offers substantially increased performance. High quality, electronically regulated membrane pumps and a new generation of valves and firmware optimizations have conferred higher precision in pressure regulation (1 mbar), lower acoustic noise, faster slew rate (~2 mBar/ms), longer lifetime and simplified calibration procedures. The pressure (300 mBar) is controlled by the PatchControl software via an USB interface.

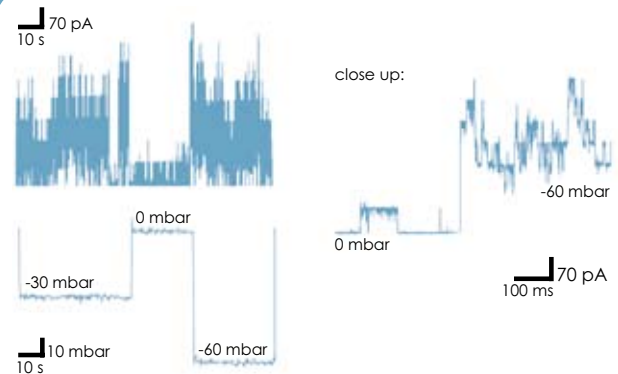
The new Suction Control Pro combines all these features with pressure clamp for accurate studies of mechano-gated ion channels, and the effect of pressure on artificial and cellular membranes.

The Suction control Pro contains an interface for communication with the analog in- and outputs of the electrophysiology recording system. This allows for precise and fast control of the applied pressure as well as the recording of the applied pressure. In this way pressure application is synchronized with voltage clamp protocols and graphically displayed together with the patch clamp data traces.

Suction Control Pro is available when purchasing a Port-a-Patch® system and is also compatible with existing Port-a-Patch® systems. Contact us if you are interested in upgrading to Suction Control Pro!



The **Suction Control Pro** offers improved performance, new experimental possibilities and is compatible with earlier versions of the Port-a-Patch® platform.



The **Port-a-Patch®** and the **Suction Control Pro** was used to study the effect of pressure on functional MscL protein reconstituted in a GUV derived planar bilayer. MscL gating was observed at a holding potential of +30mV at the indicated negative hydrostatic pressure in symmetrical recording solutions.

Europe

Nanion Technologies GmbH
Erzgiesserei-Strasse 4
80335 Munich
Germany
phone: +49 89 218997972
fax: +49 89 218997960
info@nanion.de
www.nanion.de

USA

Nanion Technologies Inc.
675 US Highway One
North Brunswick, NJ 08902
USA
phone: 1-888-9-NANION
info@naniontech.com
www.naniontech.com

nanjion