

# The Port-a-Patch<sup>®</sup> NPC<sup>®</sup>-1 Perfusion System

nanjion

Nanon Technologies GmbH



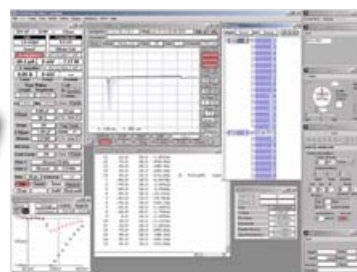
## Perfusion Chamber

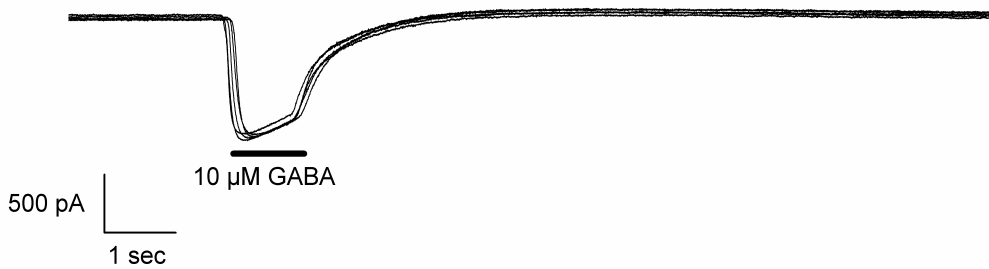
- fits all perfusion systems
- continuous flow
- easy add on



Repetitive solution exchange with the NPC<sup>®</sup>-1 perfusion chamber. High and low external potassium concentrations were applied to RBL cell expressing inward rectifying potassium channels. This figure shows that a steady state is reached in less than 200 ms.

The Port-a-Patch<sup>®</sup> is Nanion's entry level device for automated electrophysiology.





Nanon's perfusion system enables computer controlled solution changes. In the recording shown on the left 10  $\mu\text{M}$  GABA were applied for one second to HEK 293 cells expressing GABA<sub>A</sub> ( $\alpha_1, \beta_2, \gamma_2$ ) receptors. An overlay of 5 repeats is displayed.

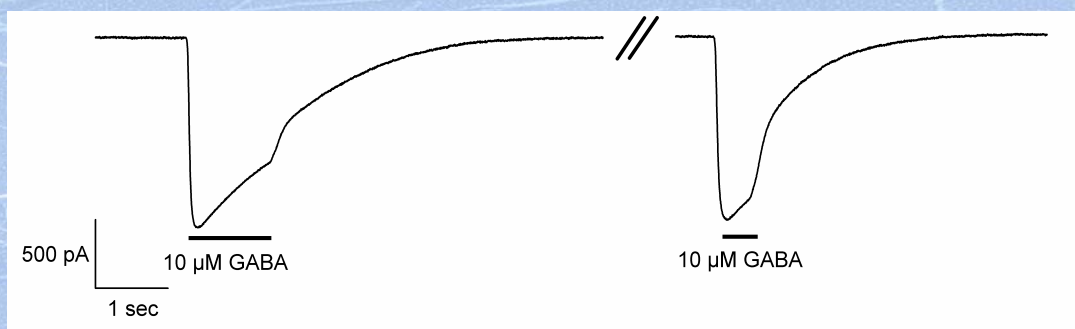


## Perfusion System

- software controlled
- stand alone solution
- suction pump included

Nanon's perfusion system provides a fast and continuous perfusion. It allows accurate application of up to eight different compounds and solutions via the patch clamp software. PatchMaster by HEKA or PatchControl 4 by Nanon can be used to control the perfusion system without an additional interface.

10  $\mu\text{M}$  GABA was applied to HEK293 cells expressing GABA<sub>A</sub> ( $\alpha_1, \beta_2, \gamma_2$ ) receptors. The first application of GABA lasted for 1.2 seconds, the second one for 500 ms. The data reveal the different time constants of desensitization and wash-out of GABA.



Smart tools for ion channel research

**Europe/USA**  
Nanon Technologies  
Erzgießereistr. 4  
80335 München  
Germany

phone: +49 89 2189 97972  
fax: +49 89 2189 97960  
info@nanion.de  
http://www.nanion.de

**Korea/China**  
Quantum Design Korea  
517-18 Kyungbin Bldg. 4F,  
Dogok-dong, Kangnam-gu,  
Seoul, Korea 135-270

phone: +82 2 2057 2710  
fax: +82 2 2057 2712  
qdk@qdkorea.com/info@qdchina.com  
www.qdkorea.com/www.qd-china.com

**Japan**  
Bio Research Center Co., Ltd.  
2-28-24, Izumi, Higashi-ku  
Nagoya 461-0001  
Japan

phone: +81 52 932 6421  
fax: +81 52 932 6755  
ohtsukik@nanion.de  
http://www.brck.co.jp

**India**  
Medi Analytika India Pvt. Ltd  
6, Adyar Bridge Road  
Adyar, Madras - 600 020  
India

phone: +91 44 2446 0988  
fax: +91 44 2446 3931  
mediana@vsnl.com  
http://www.medianalytika.com

**Taiwan**  
Sunpoint Scientific Instrument  
9f-1, No.146, Wen Shing Rd.  
Taoyuan Hsien 333  
Taiwan, R.O.C

phone: +886 3 327 3889  
fax: +886 3 327 3908  
sunp55@ms36.hinet.net  
http://www.sun-point.com.tw