



The **Port-a-Patch[®]**

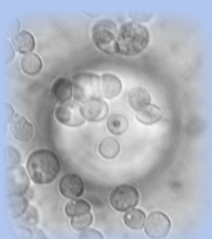
1
easy to use
complete
patch clamp
setup

2
planar patch
clamp chips

3
fast fluid
exchange

4
small compound
consumption

5
high quality
recordings



PatchControl 4.0 now available:

Get a closer look at the features of new PatchControl 4.0



The Port-a-Patch[®] has been successfully introduced as a high-quality automated patch clamp platform for industry and academia.

Nanion is now launching the new version of the Port-a-Patch[®] software: *PatchControl 4.0*. In addition to controlling Gigaseal formation, the patch clamp experiment can now be completely pre-programmed.

New features



More automation. Powerful macros for enhanced automation of experiments. Designing experiments is quick and simple.



More visualisation. The workflow window in the log file offers more space to add information describing the things you need.



More users. Multi-user functionality enabling every user to create their own logfile and macros.



More interface. The graphical user interface is more user-friendly.



More compatible. Windows Vista compatible. *PClamp 9* and *10*; *Pulse* and *PatchMaster* compatible. *PatchControl 4.0* is now compatible with more operating systems, more electro-physiology and most importantly...you!



The
Port-a-Patch®

1
easy to use
complete
patch clamp
setup

2
planar patch
clamp chips

3
fast fluid
exchange

4
small compound
consumption

5
high quality
recordings

As an example of one of the improved functionalities, new macros have been designed to incorporate automated internal perfusion control.

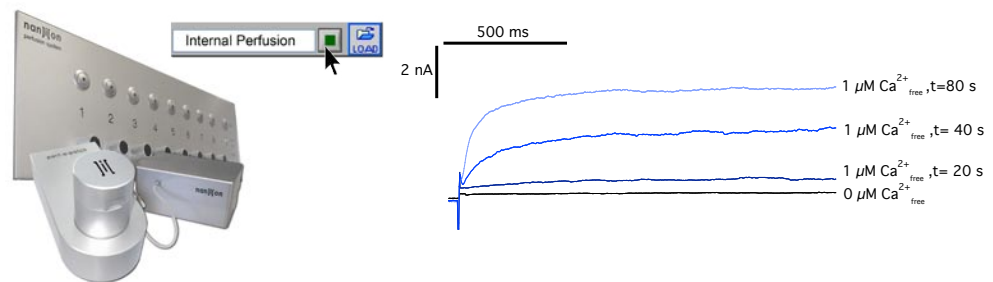


Fig. 1 **Whole cell recordings performed with the Port-a-Patch® using automated internal perfusion.** Currents mediated by BK channels (large conductance Ca^{2+} and voltage-activated K^+ channels) expressed in CHO cells. Currents were elicited by a voltage-step from -80 mV to $+80$ mV before ($0 \mu\text{M Ca}^{2+}$); and 20, 40 and 80 s after the internal application of $1 \mu\text{M free Ca}^{2+}$.

New *PatchControl 4.0* will be released at the Neuroscience conference in San Diego, CA, November 3 - 7 2007. Please visit us at booth no. 613. All Port-a-Patch users are welcome to take a CD containing *PatchControl 4.0*.

For more information or if you have any questions please see www.nanion.de or contact us at info@nanion.de.