

# PatchControl 384/DataControl 384

## Version 1.6: High throughput. Made easy.

- New version of PatchControl 384 & DataControl 384
- Two modes of action for the Biomek - standard and advanced
- New online analysis functions
- Simplified quality control features
- Easy experimental set-up
- Scanning and storage of NPC-384 barcode
- New display features

### PatchControl 384/ DataControl 384 Version 1.6

#### Improving usability

The new release of the SyncroPatch software includes an update of both PatchControl 384 and DataControl 384 for efficient recording and analysis of data recorded on the SyncroPatch 384/768PE. The keyword for this software release is **simplicity**. There are now two modes of action for controlling the Biomek - 'standard' and 'advanced'. The standard mode has been designed with simplicity in mind, minimizing the number of commands required to set up an experiment. Creating a method from scratch has also been made a lot easier. In advanced mode, the same method can be edited with additional commands, thus maintaining flexibility. New recording and analysis features include '**Ligand Puff**' command for more robust and reproducible recordings of ligand-gated ion channels; **exponential fits of raw traces** for estimation of kinetics; **reversal potential**; **slope and decay analysis**, amongst others.



The SyncroPatch 384PE

#### Key improvements of the PatchControl 384 software

- Two modes of action: Standard and advanced
- Clearer user interface - Simplified experimental set-up
- New 'Ligand Puff' command to minimize desensitization of ligand-gated ion channels
- New online analysis functions: reversal potential, time to peak, tau, slope and more
- New display features, e.g. select wells in a row

PatchControl 384  
DataControl 384

nanjion

**Nanion Europe**

info@nanion.de

Phone: +49 89 2190 95-0

www.nanion.de

**Nanion USA**

info@naniontech.com

Phone: 1-888-9-NANION

**Nanion China**

info@nanion.cn

phone: +86 1082 17 63 86

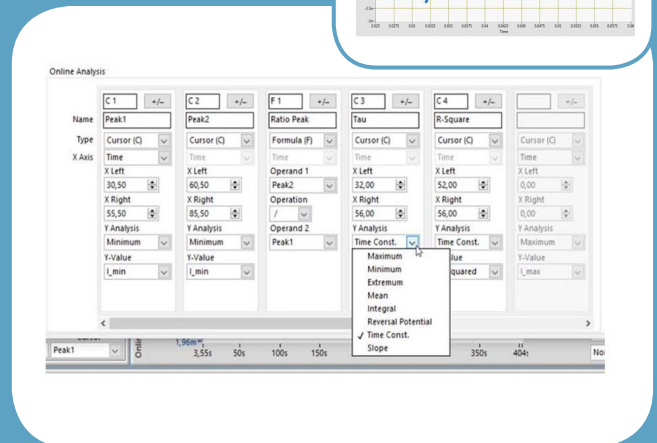
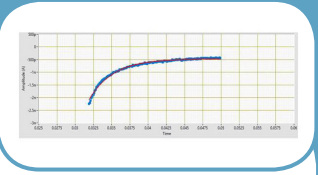
**Nanion Japan**

ohtsuki@nanion.jp

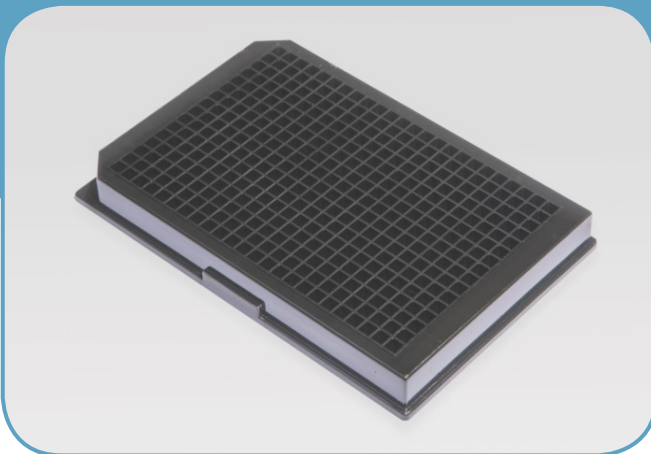
phone: +81 3 6457 8776

**nanjion**Key improvements of the **DataControl 384** software

- Revision of analysis and normalization parameters
- New online analysis functions: reversal potential, time to peak, advanced integral, tau & slope
- Display averaged data, including average online analysis and traces for easier overview of compound effects
- Improved PDF reporting
- QC summary now possible without inclusion of customer information for easier support and troubleshooting



New online analysis functions, e.g. exponential fit for tau calculation



Every NPC-384 chip has a barcode and this is scanned at the beginning of every experiment for easier tracking and support.

This upgrade is available to all SyncroPatch 384PE customers **free-of-charge** with an annual service contract. Want to find out more about how to upgrade? Please contact: [support.384PE@nanion.de](mailto:support.384PE@nanion.de) for more information.

## PatchControl 384/DataControl 384: version 1.6 update

*“The improved current trace colour co-ordination in the PDF reports as well as the new feature which shows the average current time-course and the average current traces helps to quickly create a clear picture of the overall test compound effect without the need to look at each individual cell. The simplified assay set-up software is much more user friendly, which has improved our experimental efficiency as it has allowed us to adjust each parameter for our wide range of cell lines and assays with greater ease.”*

SB Drug Discovery Electrophysiology Group

## PatchControl 384/DataControl 384: Version 1.6.