

DataControl 384

SyncroPatch 384

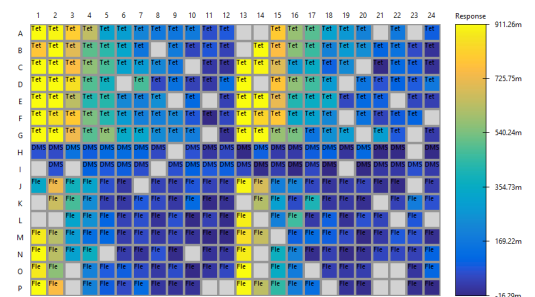
Effortless analysis with DataControl 384

DataControl 384 is a standalone program designed for swift and effortless loading and analysis of data recorded on the SyncroPatch 384.

Key benefits:

- **Online Analysis:** Monitor and analyze relevant parameters during the recording.
- **Automated IV analysis:** Automatic IV construction and determination of half-maximal activation or inactivation potentials (V_{half}).
- **Analysis of compound effects:** Automated concentration response curve generation and IC_{50}/EC_{50} determination.
- **Control Well Correction:** Automatically correct for rundown for more accurate IC_{50} and EC_{50} values.
- **Data Report:** Export raw data or summary of results and experimental details in table, word or pdf format. Export graphs as jpg, tif, png, eps, pdf or svg files.

Flexible and powerful analysis at your fingertips



Heat maps:

- Observe distribution of V_{half} or compound effects.
- Easy color-coded overview.
- High resolution export of heat map view.

Contact us today

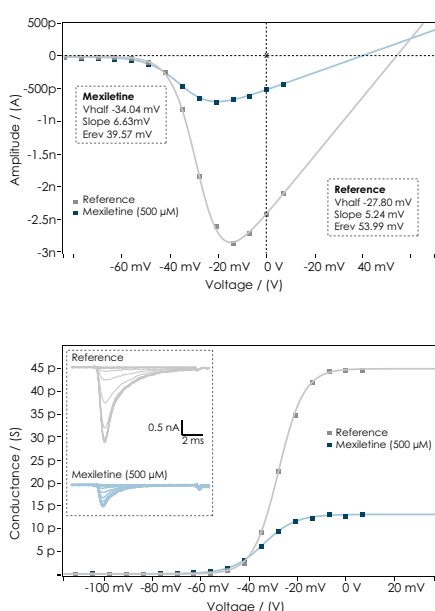


DataControl 384 for the SyncroPatch 384

Analyze all your data with a few clicks

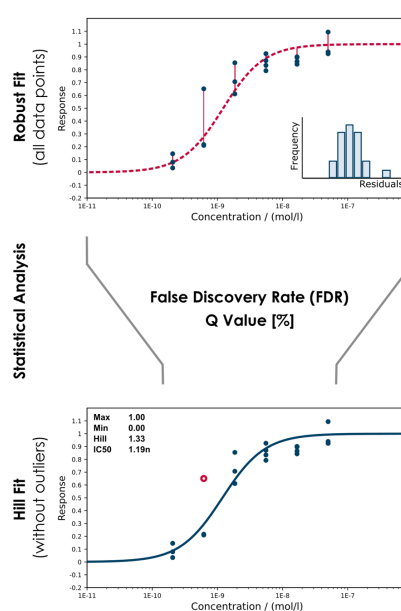
Automated IV analysis

- Individual and averaged IV curves: 3 modes of automated IV fitting (individual, averaged and pooled).
- Precise fit with automated quality control.
- Generate accurate activation and inactivation curves pre- and post-compound.
- Automatically measure effects of state-dependent compounds on V_{half} and G_{max} .
- Optimized for all voltage-gated ion channels (activation, inactivation, tail currents).



CRC analysis with outlier detection

- Use pre-defined templates to construct concentration response curves.
- Automatically determine IC_{50} values of inhibitors or EC_{50} values of positive modulators (activators, potentiators).
- Automatically detect and eliminate outlier points for improved curve fitting.
- Automated QC filters out unsuccessful wells based on user defined parameters for faster analysis.



Control well correction

- Use vehicle control wells to determine rundown effect over time.
- Automatically subtract rundown from compound wells for more accurate calculation of concentration response.
- Turn on and off control well correction with a simple click.

