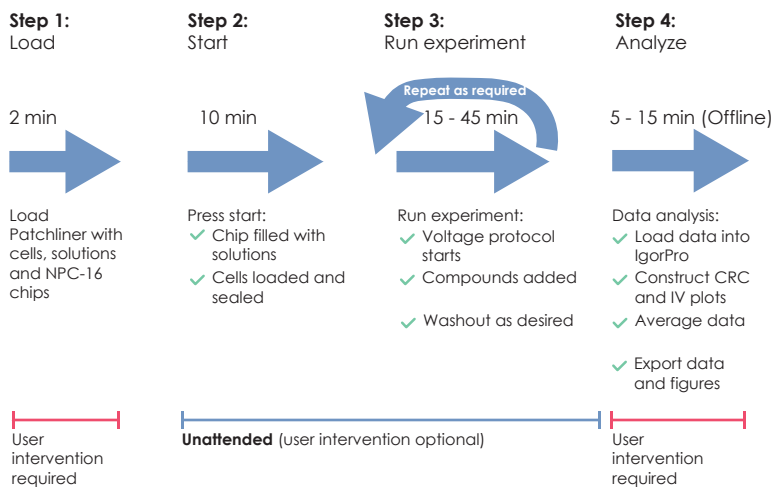


PatchControlHT Suite - Patchliner



Complete software suite for the Patchliner

The PatchControlHT suite is a fully integrated software solution for experiment execution, data acquisition and data analysis on the Patchliner.



Key benefits:

- **Full automation of experiment execution** regardless of experimental design.
- **Retain control** for changes on the fly.
- **User and Expert modes of operation** for running routine experiments or designing new and sophisticated protocols.
- **PatchMaster software** for data acquisition.
- **Automated analysis in Igor Pro:** for concentration response curves and IC/EC₅₀ calculations, as well as IV curves with V_{half} of activation/inactivation values.



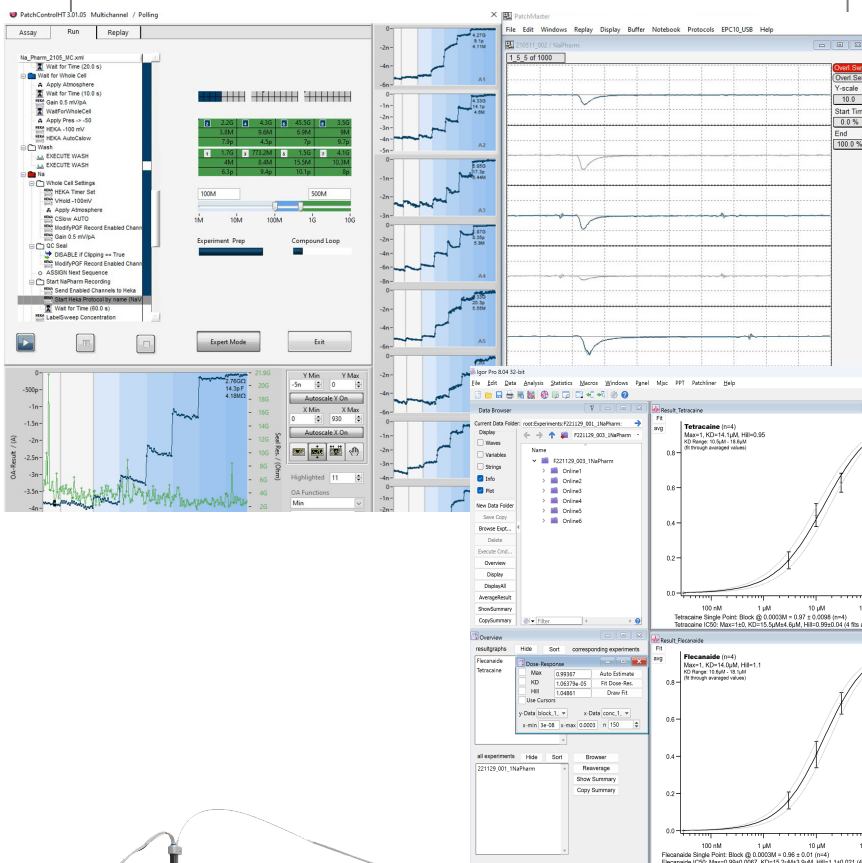
Contact us today

Software suite for the Patchliner

Complete data acquisition and analysis

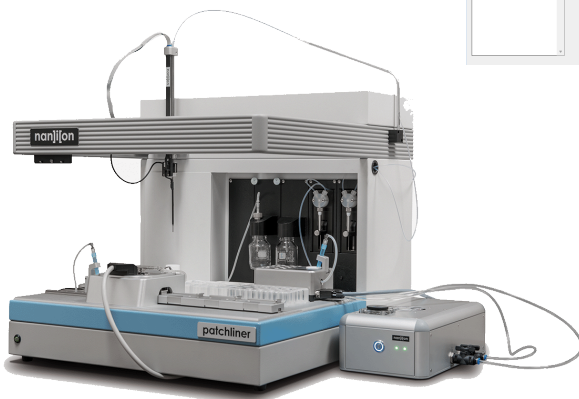
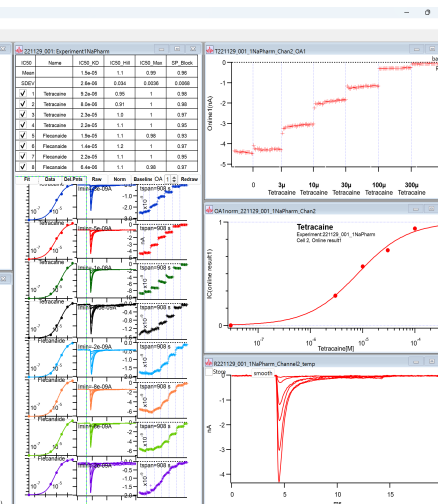
PatchControlHT

- Simply drag and drop robot and recording commands into the list.
- Commands logically grouped in folders for easy overview
- Fully automated procedures.
- Changes on the fly are possible.
- Color-coded overview of cell parameters and timecourse of experiment.



PatchMaster

- HEKA EPC-10 amplifier(s).
- 4 or 8 amplifier channels.
- Whole cell or on cell mode.
- Voltage clamp, current clamp, dynamic clamp
- Raw current traces displayed.
- Easy transfer of pulse protocols and online analysis settings from your manual rig.



Igor Pro

- Easy import of data.
- Automated construction of IVs and CRCs.
- Pool data from multiple NPC-16 chips.
- Export of graphs and data in different formats.