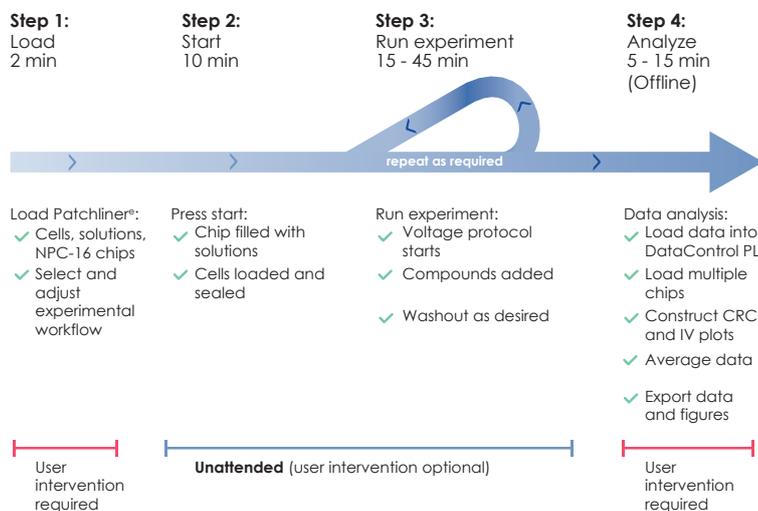


PatchlinerControl Patchliner^e

Simple, comprehensive experiment setup and data acquisition

PatchlinerControl is a proprietary software for setup of experiments and data acquisition for the Patchliner^e. Fully automated workflows, QC, and real-time data analysis ensure maximum efficiency.



Key benefits:

- **Fully automated experiment execution.**
- **Real-time control** and on-the-fly adjustments.
- **Fast, intuitive setup** with predefined commands.
- **Clear, live data visualization** of individual wells.
- **Automated quality control** for reliable data acquisition and analysis.
- **Advanced analysis in DataControl PL:** CRCs, IC₅₀/EC₅₀ and IV curves with V_{1/2} values.

PatchlinerControl

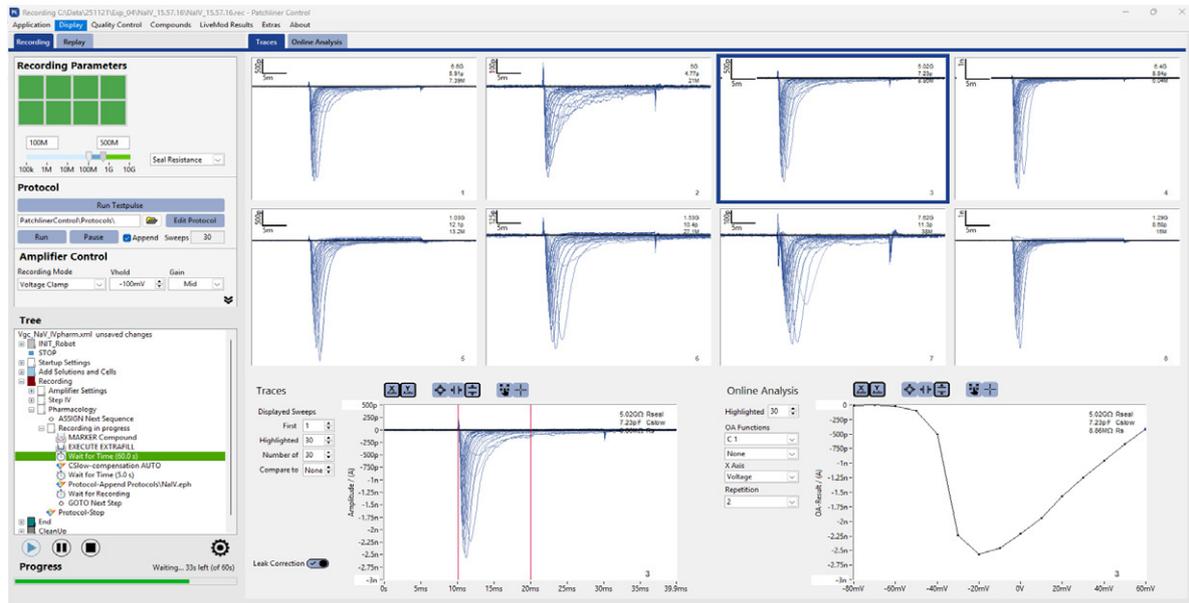
Easy data acquisition and online analysis

Chip and experiment overview

- Color-coded chip map for quick visualization of success rates.
- Live recording mode plus replay of recorded data.
- Load, edit, and run voltage protocols with ease.

Real-time data overview

- Display active recording or pre-recorded data.
- View all 8 wells at a glance.
- Switch instantly between raw traces and online analysis.



Experiment setup and execution

- Choose from a library of commands for simple yet powerful experiment setup.
- Track experiment progress at a glance at any timepoint.
- Stop, pause, or restart the experiment anytime.
- Make changes on-the-fly during execution.
- Seamless transfer of voltage protocols across Nanion APC platforms.

Quick online analysis

- View real-time analysis for any individual well.
- Select and display the most relevant analysis parameters.

Contact us today

