

# PATCHLINER

## PatchControl HT

nanjion

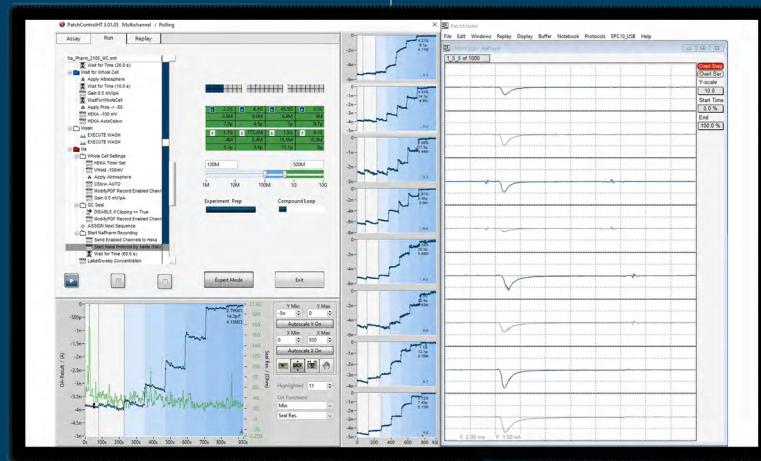
### Advanced online analysis display

- Solution additions labeled in online analysis graphs
- Switching between different online analysis parameters during recording
- Zoom-in graph of a selected cell
- Interface adapted to SyncroPatch software for easy transition between systems
- Easy online monitoring and post-hoc review of experiment

### PatchControl HT

#### Simplified user mode interface in addition to expert mode

- One-click switch between expert mode and user mode (and vice versa)
- Easy operation via tabs for assay selection, running an experiment and replay of data
- Minimal training required for operating the Patchliner
- Streamlined overview of experimental parameters
- Adjustable color scale for seal resistance display
- Full flexibility for sophisticated assay development



Watch me in action



[www.nanion.de](http://www.nanion.de)

The Patchliner is a fully automated planar patch clamp instrument recording from up to 8 cells simultaneously. With its vast experimental freedom and gigaseal data quality, the Patchliner is one of the most versatile patch clamp instruments on the market.



Patchliner with Dynamite<sup>8</sup>



Nanion Technologies GmbH  
Ganghoferstraße 70a  
80339 München / Munich

Telefon: +49 / (0)89 2190 95 - 072  
Telefax: +49 / (0)89 2189 97 - 960  
E-Mail: [info@nanion.de](mailto:info@nanion.de)