

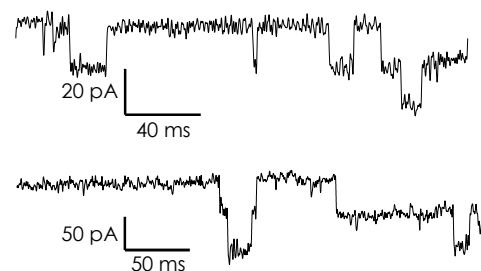
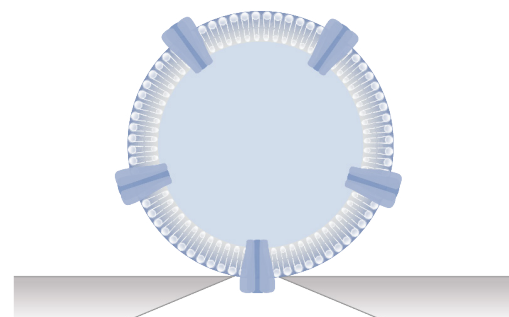


Single Channel Recordings Patchliner

Biophysical properties of single channels on the Patchliner

Automated patch clamp devices are typically used for routine screening of compounds and target validation, but what if you want to get down to the nitty gritty biophysical properties of single channels? With the Patchliner, a fully automated patch clamp system for recording up to 8 cells simultaneously, you can do just that! The Patchliner uses high quality HEKA EPC10 amplifiers, often used in manual patch clamp rigs, coupled with borosilicate glass chips for low noise, low capacitance recordings from up to 8 cells at a time. Use NPC-16 high resistance chips and control suction parameters manually to make sure you get high giga-Ohm seals in on-cell mode for single channel recordings.

Want to do whole cell recordings after all? No problem, simply use suction to break into the whole cell mode and continue your recording.



Contact us today

Patchliner

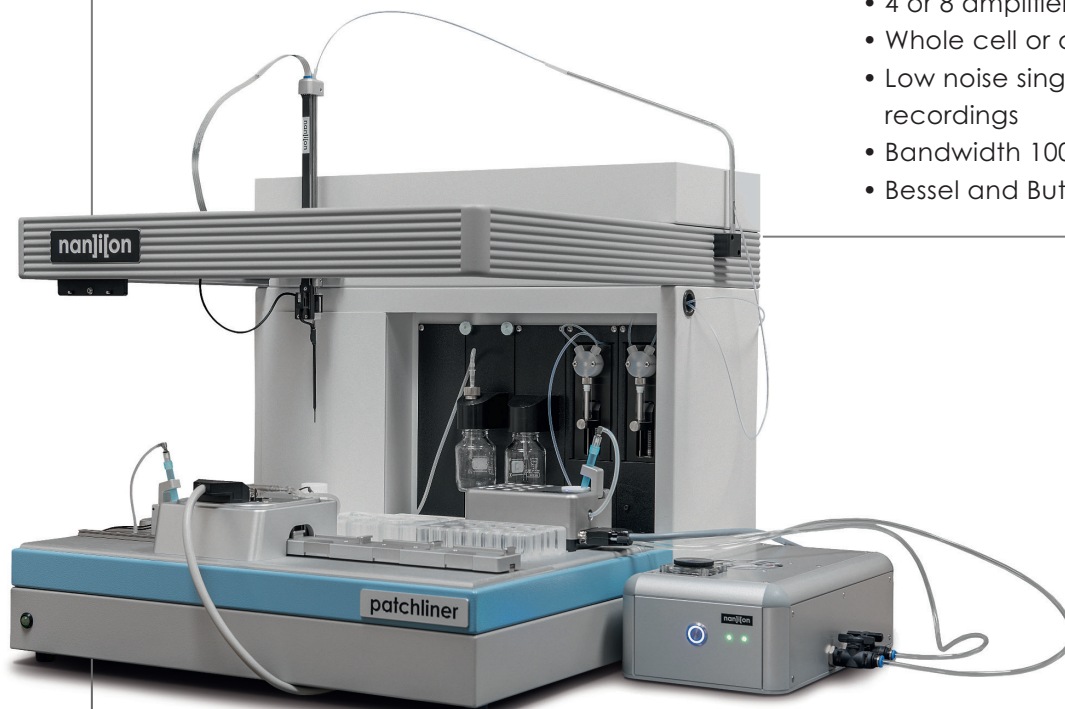
Multiple recording configurations

Borosilicate glass chips

- Low capacitance
- Low noise
- Long-lasting recordings
- High resistance for cell attached recordings

Amplifier

- HEKA EPC-10 amplifier(s)
- 4 or 8 amplifier channels
- Whole cell or on cell mode
- Low noise single channel recordings
- Bandwidth 100 kHz
- Bessel and Butterworth filters



Applications

- Voltage-, ligand- and heat-activated ion channels
- Mechanical stimulation using shear stress
- Pressure activation using suction
- Single channel recordings in cell-attached mode
- Automated or manual control of suction