



Axon[™] Patch-Clamp

The Axon Instruments™ portfolio provides comprehensive solutions for patch-clamping that includes amplifiers, digitizer, software, and accessories. Our best-in-class instruments facilitate the entire range of patch-clamp technique experiments from the smallest single channel to the largest macroscopic recordings. The addition of Axon pCLAMP™ 11 Software Suite creates a streamlined workflow, allowing for sophisticated and efficient experiments, and higher quality data generation.

Axon Instruments



Acquisition—get more flexibility and easier acquisition

- Enhanced protocol editor in pCLAMP 11 Software allows for longer and more sophisticated protocols to be programmed and reduces protocol setup time
- · Multiple windows can be simultaneously displayed during a Membrane Test protocol, making it easier to patch multiple cells
- Gap-free mode allows the programming of command and digital outputs in a continuous recording, allowing you to enhance
 protocol design and trigger actions within the protocol automatically



Amplification—amplify your success

- · Control via pCLAMP software maximizes flexibility in setting up and performing different experimental protocols
- · A wide range of amplifiers means we can offer the ideal choice for your application, providing the best possible data
- Our lowest-noise amplifier/digitizer instruments provide the highest signal resolution, and eliminate 50 or 60 Hz line-frequency noise contamination



Digitization—get better data, faster

- High rate of sampling (500kHz) means a more faithful digitization process, allowing for precise digital reconstruction of the analogue signal due to minimal information loss
- · Eight channels can be sampled and digitized simultaneously, increasing data acquisition speed
- The HumSilencer[™] option with the Digidata range eliminates of 50 and 60Hz line-frequency noise, making the biological signal clearer without distorting the data



Analysis—discover more while saving time

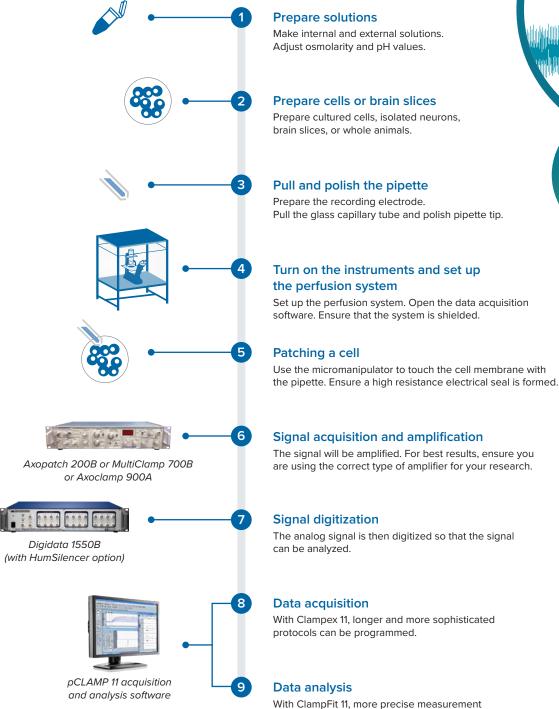
- Acquisition and analysis are combined in one package, streamlining the entire process
- Advanced modules allow action potential and population spike measurements to be made automatically
- New Batch Analysis feature allows multiple datasets to be analysed using an identical macro; saving time on data analysis and ensuring the exact same analysis is carried out on each dataset

pCLAMP is the most cited patch-clamp software!

Which amplifier is right for me?

	Axopatch™ 200B Amplifier*	MultiClamp™ 700B Amplifier	Axoclamp™ 900A Amplifier
Single-channel recording	• • •	•	
Whole-cell voltage-clamp	• • •	• • •	
Whole-cell current-clamp	•	• • •	• • •
Bilayer study	• • •	• • •	
Extracellular field-potential recording	•	• • •	• • •
Amperometry/voltammetry study	• •	• • •	
Nanopore study	• • •	•	
Intracellular sharp-electrode recording	•	• • •	• • •
Two-electrode voltage-clamp recording			• • •

Electrophysiology workflow using **Axon Instruments**



and faster data analysis are now possible.

with HumSilencer™

Raw Data

HumSilencer
eliminates
50/60 Hz
line-frequency

noise

Fast and effective data analysis turns data into actionable insights and publication-ready results.

Applications

Our solutions can be used across almost the entire range of electrophysiology techniques, from single channel to whole cell to extracellular field-potential recording.



Extracellular recording

- · Single-unit recording
- · Multi-unit recording
- Field-potential recordings
- · Amperometry/voltammetry



Intracellular recording

- · Cell-attached/excised patch-clamp
- Whole-cell voltage-clamp
- · Whole-cell current-clamp
- · Sharp-electrode



University of Texas

uses Axon to evaluate signaling mechanisms in the brain underlying memory and recall.

Allegheny College uses Axon to investigate how amyloid beta peptide block ion channels in Alzheimer's disease.





University of Michigan uses Axon to investigate NMDAR

receptor blockers.



Want to learn more about electrophysiology techniques? Download the free Axon Guide.



Contact Us

Phone: +1.800.635.5577

Web: www.moleculardevices.com

Email: info@moldev.com

Check our website for a current

listing of worldwide distributors.

Regional Offices

 USA and Canada
 +1.800.635.5577
 Taiwan/Hong Kong
 +886.2.2656.7585

 United Kingdom
 +44.118.944.8000
 Japan
 +81.3.6362.9109

 Europe*
 00800.665.32860
 South Korea
 +82.2.3471.9531

 China
 +86.4008203586
 India
 +91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Iralay, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom