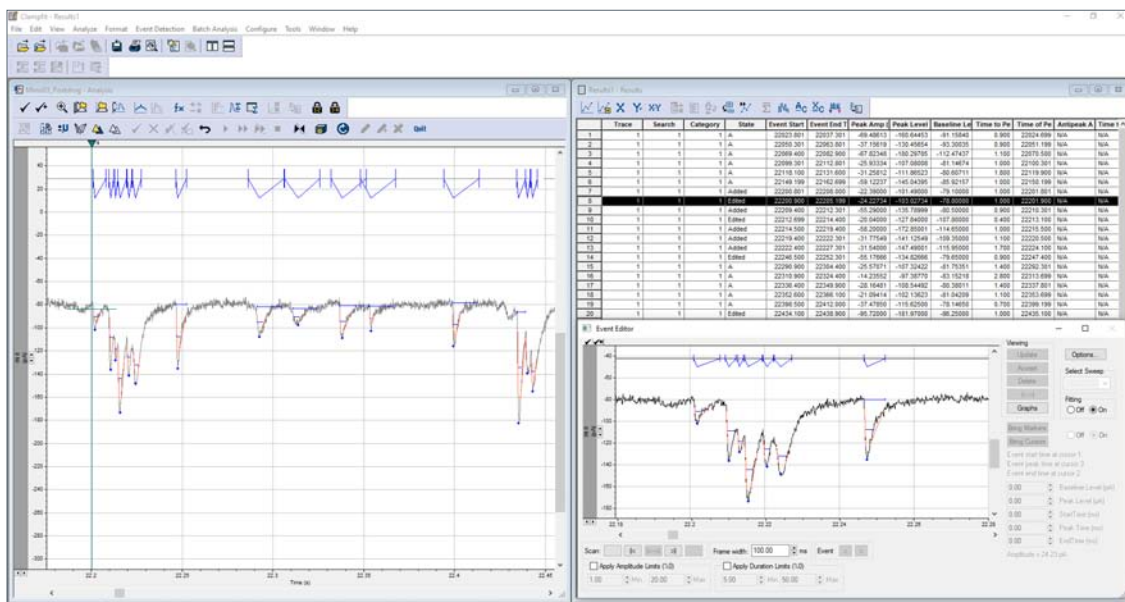


New Minis Analysis in Axon™ pCLAMP™ Software: Minis Search

The Minis Search is a new event detection feature in the Axon pCLAMP 11 software. This search is particularly designed to analyze spontaneous or miniature excitatory postsynaptic currents (sEPSCs or mEPSCs) and spontaneous or miniature inhibitory postsynaptic currents (sIPSCs or mIPSCs). It provides a streamlined workflow for detecting minis events acquired in a continuous or episodic-sweep mode. In addition, it enables accurate measurements of detected events and simplifies the process of event addition and deletion. Overall, this comprehensive analysis allows neuroscientists to analyze minis events in an easier, faster, and more accurate way.



Key functions

- Streamline the workflow for detecting minis events
- Enable accurate measurement of each detected event
- Easy to add a missing event
- Easy to delete an accepted event
- Able to detect superimposed events
- Able to adjust the event baseline

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
United Kingdom +44.118.944.8000
Europe* 00800.665.32860
China +86.4008203586
Taiwan/Hong Kong +886.2.2656.7585
Japan +81.3.6362.9109
South Korea +82.2.3471.9531
India +91.73.8661.1198

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

