



# SoftPanel Controller

## KEY FEATURES

- Offers a more conventional interface to Axon Multiclamp 700B and Axoclamp 900A microelectrode amplifier controls
- Physically replicates the essential functions of the Commander Software
- Predefined color-coded magnetic overlay for both voltage clamp and current clamp

The SoftPanel Controller provides a more conventional interface to amplifier control. Designed as a hardware extension to replicate essential amplifier function for Axoclamp™ and MultiClamp™ Commander Software, the SoftPanel Controller includes 8 configurable knobs and 8 configurable buttons. Using these knobs replicate continuous mouse controls (“gliders”), while the buttons replicate single-click mouse controls without negating the many benefits afforded by computer control of the amplifier. The SoftPanel Controller can be easily set up with the computer via USB 2.0 connection.

## Specifications

Dimensions (in.)	3.7 (H) x 5 (W) x 7 (L)
Dimensions (cm)	9.5 (H) x 12.7 (W) x 17.8 (L)
Weight (lbs.)	1.70
Compatibility	Axoclamp 900A and Multiclamp 700B Amplifiers
Communications and power	USB-2 port compatible cable with USB type A to USB type B connections (provided)
Includes overlays for keys and buttons	For both Axoclamp 900A or Multiclamp 700B Amplifiers



The Softpanel Overlay for the MultiClamp 700B Microelectrode Amplifier.

The Softpanel Overlay for the Axoclamp 900A Microelectrode Amplifier.

**Contact Us**

Phone: +1-800-635-5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto:info@moldev.com)  
 Check our website for a current listing of worldwide distributors.