

PowerPac™ Adaptor Catalog Numbers 164-5062, 164-5064



Warning:

Bio-Rad PowerPac power supplies and electrophoresis cells are designed and certified to conform to IEC 61010-1, an internationally accepted standard for electrical safety of laboratory instrumentation. Compliant products are safe to use when operated in accordance with manufacturer's instructions. The safety certifications of our power supplies do not, however, extend to electrophoresis cells or other accessories connected to them.

This PowerPac adaptor is designed to IEC 61010-1 specifications for connection isolation and flammability. It is provided to allow customers to use our power supplies with some other manufacturer's cells and accessories. Note, however, that use of this adaptor voids our PowerPac power supply IEC 61010-1 safety certification and may present a safety hazard to the user depending on the safety of the equipment being attached. Please check with the electrophoresis cell or other accessory manufacturer for compliance issues.

Instructions:

The following guidelines should be observed and followed when using the Bio-Rad PowerPac adaptor:

- Keep the PowerPac adaptor connected to the electrophoresis cell leads rather than the PowerPac power supply.
- Follow all safety precautions as listed in the safety section of the PowerPac power supply instruction manual.

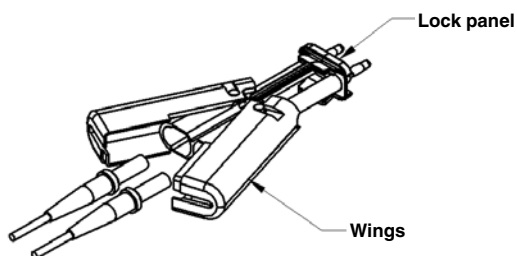


Fig. 1.

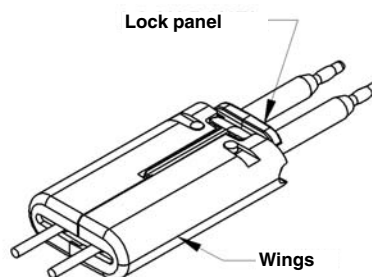


Fig. 2.

1. Slide the lock panel up to open the PowerPac adaptor wings. See Figure 1.
2. Insert the leads for the non certified electrophoresis cell into the PowerPac adaptor. Ensure that the polarity of the cables is maintained (i.e., red to red and black to black).
3. Close the PowerPac adaptor wings and slide the lock panel down. See Figure 2.
4. Plug the PowerPac adaptor into the correct output terminals of the PowerPac power supply maintaining the polarity of the cables (i.e., red to red and black to black).