

## Procedure to Remove the Power Cable Tie-Bar

**Note:** This removal voids IEC1010-1\* electrical safety standards.

These Bio-Rad electrophoresis power cables are joined together by a tie-bar that insures compliance to the international IEC1010-1 electrical safety standard. However, the tie-bar on the cables can be removed in order to fit power supplies that do not conform to IEC1010-1 regulations. The tie-bar is removed by cutting the four retention tabs from the cable leads and sliding off the bar. Removal of the retention tabs is permanent and voids the IEC1010-1 safety certification.

\*IEC1010-1 is an internationally accepted electrical safety standard for laboratory instruments.