

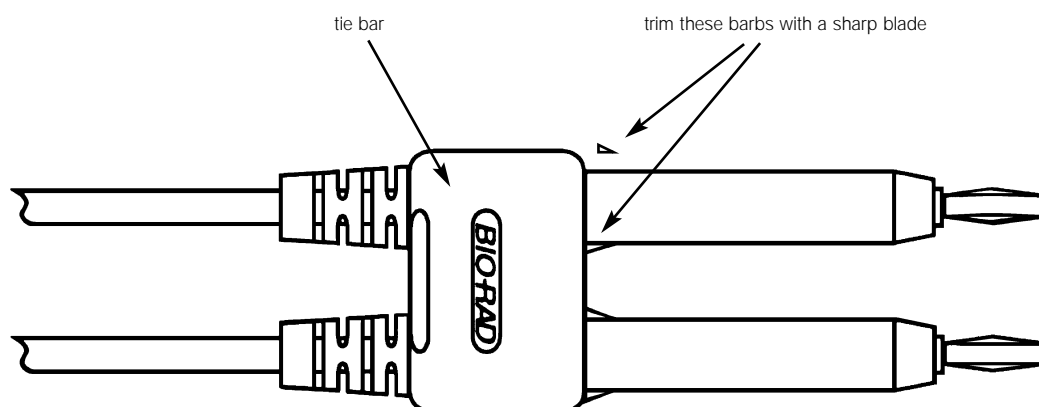
Removable Tie Bar on EN61010-1 Duplex Power Cables

Bio-Rad Laboratories, Life Science Group, Technical Service,
2000 Alfred Nobel Drive, Hercules, CA 94547

The tie bar on the most recent version of EN61010-1 duplex power cables can be removed (see figure below). This option is only possible with cables whose plugs are not fixed into the tie bar, as indicated by the ability to rotate the plugs by twisting the leads.

The tie bar can be removed from the cables by trimming the barbs with a razor blade or scalpel.

Removing the tie bar is not recommended by Bio-Rad and will void the EN61010-0 certification. This option is available because some power supplies will not work with Bio-Rad's EN61010 certified connector.

**BIO-RAD****Bio-Rad
Laboratories**

Life Science
Group

Web Site www.bio-rad.com U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
China 86-10-62051850/51 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France 01 43 90 46 90 Germany 089 318 84-0
Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 39-02-216091 Japan 03-5811-6270 Korea 82-2-3473-4460
Latin America 305-894-5950 Mexico 514-2210 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 22-74-18-70 Russia 7 095 979 98 00
Singapore 65-2729877 Spain 34-91-661-7085 Sweden 46 (0)8-55 51 27 00 Switzerland 01-809 55 55 United Kingdom 0800-181134