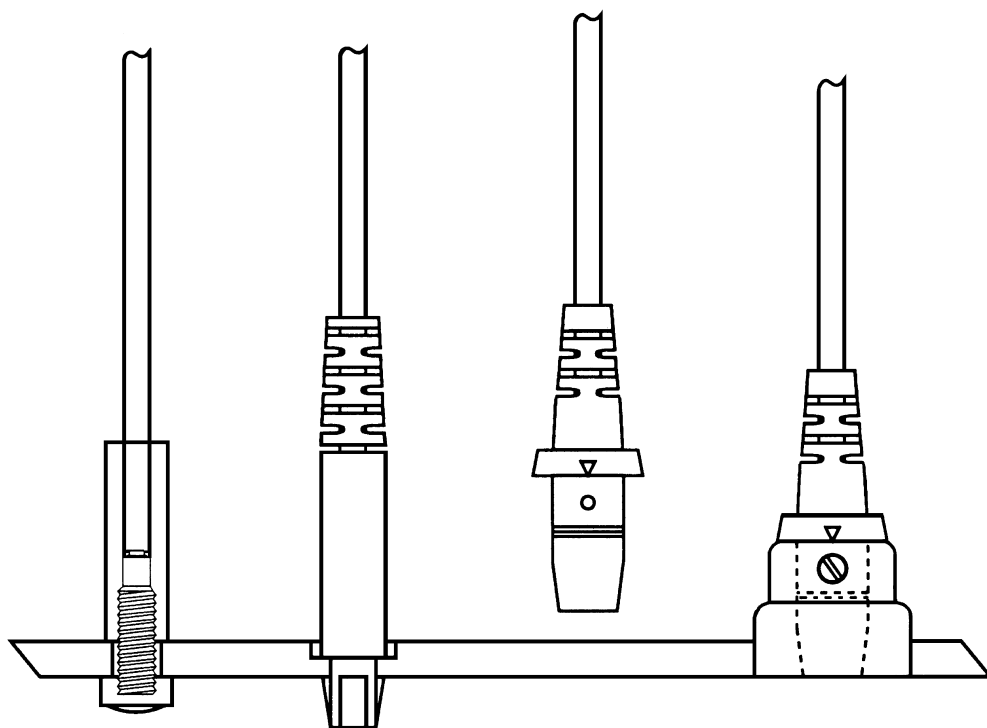


Electrophoresis Cable Plug Ends

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

Three different types of electrophoresis cable plug ends are available from Bio-Rad (see figure below). The old-style screw-on jack and the molded plastic plug ends work with 21 mm banana plugs. The EN61010 plug end is for longer 37 mm banana plugs.



Old-style screw-on jack
catalog #165-1805

Molded plug end in
recessed hole
catalog #165-1915
catalog #165-2947

EN61010 plug end
catalog #165-2948

EN61010 plug end
in socket on lid

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