

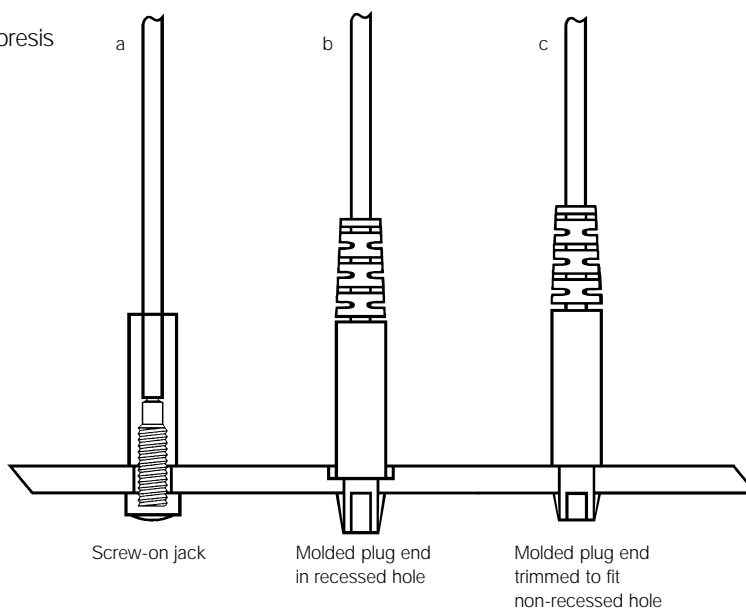
Modification of Molded Plug End to Fit Lids with Non-Recessed Holes

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

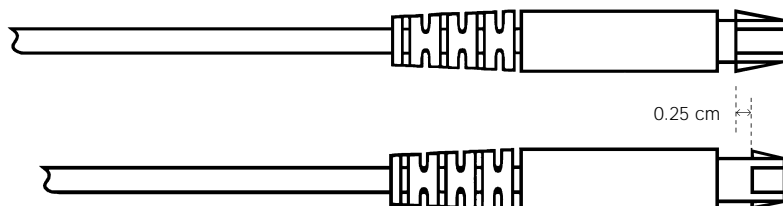
The molded plug end on replacement electrophoresis cables (catalog #165-2947, straight wires, or catalog #165-1915, coiled wires) can be easily modified to fit electrophoresis lids without recessed holes (see figure at right).

For comparison purposes, the screw-on jack (a) and the molded plug end (b) are shown.

A molded plug end that has been trimmed to fit a non-recessed hole is also shown (c).



To modify a molded plug end to fit a lid with non-recessed holes, trim all four barbs on the molded plug end with a razor blade or scalpel, leaving a space between the barbs and the base of the plug end equal to the thickness of the lid. Typically, trimming about 0.25 cm from the barb ends is sufficient (see figure below).



Trim all four barbs on the plug end with a sharp blade

To order replacement cables, phone 1-800-4BIORAD within the US, or contact your local Bio-Rad representative.



**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