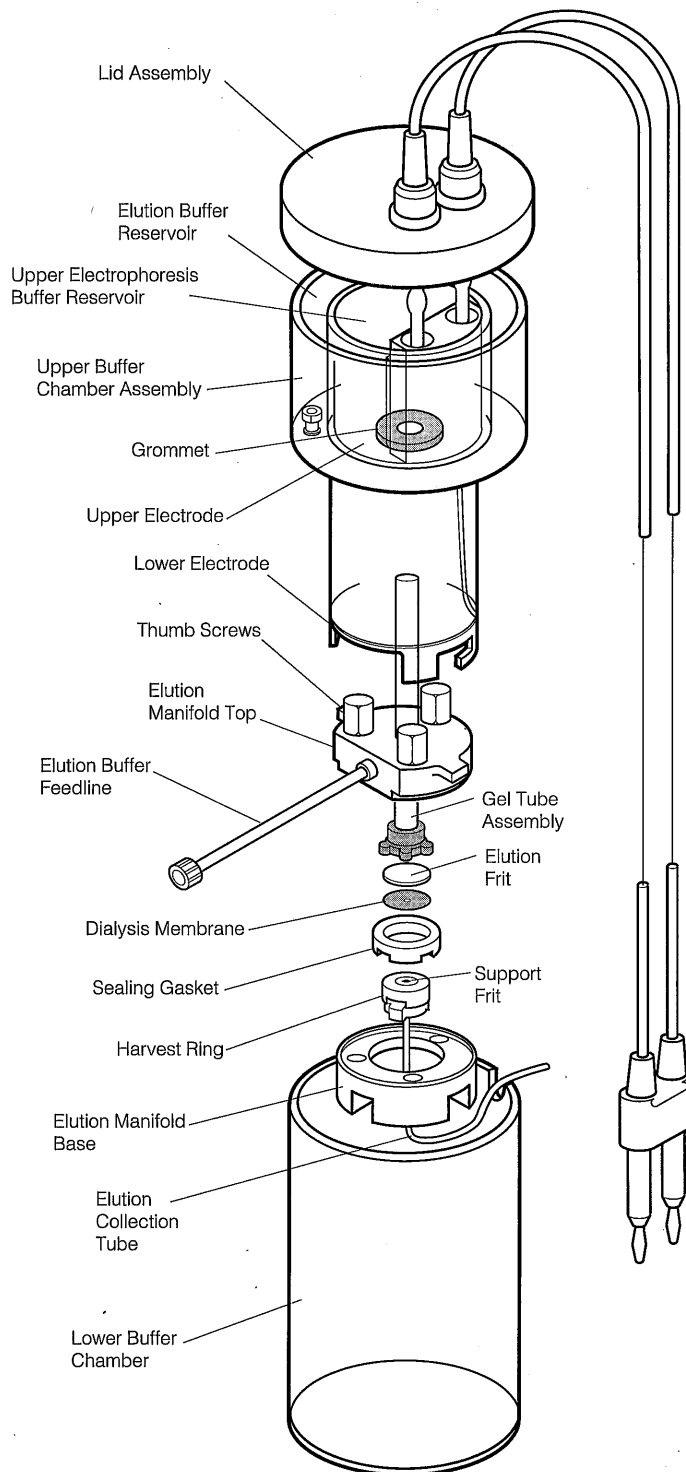




Mini Prep Cell

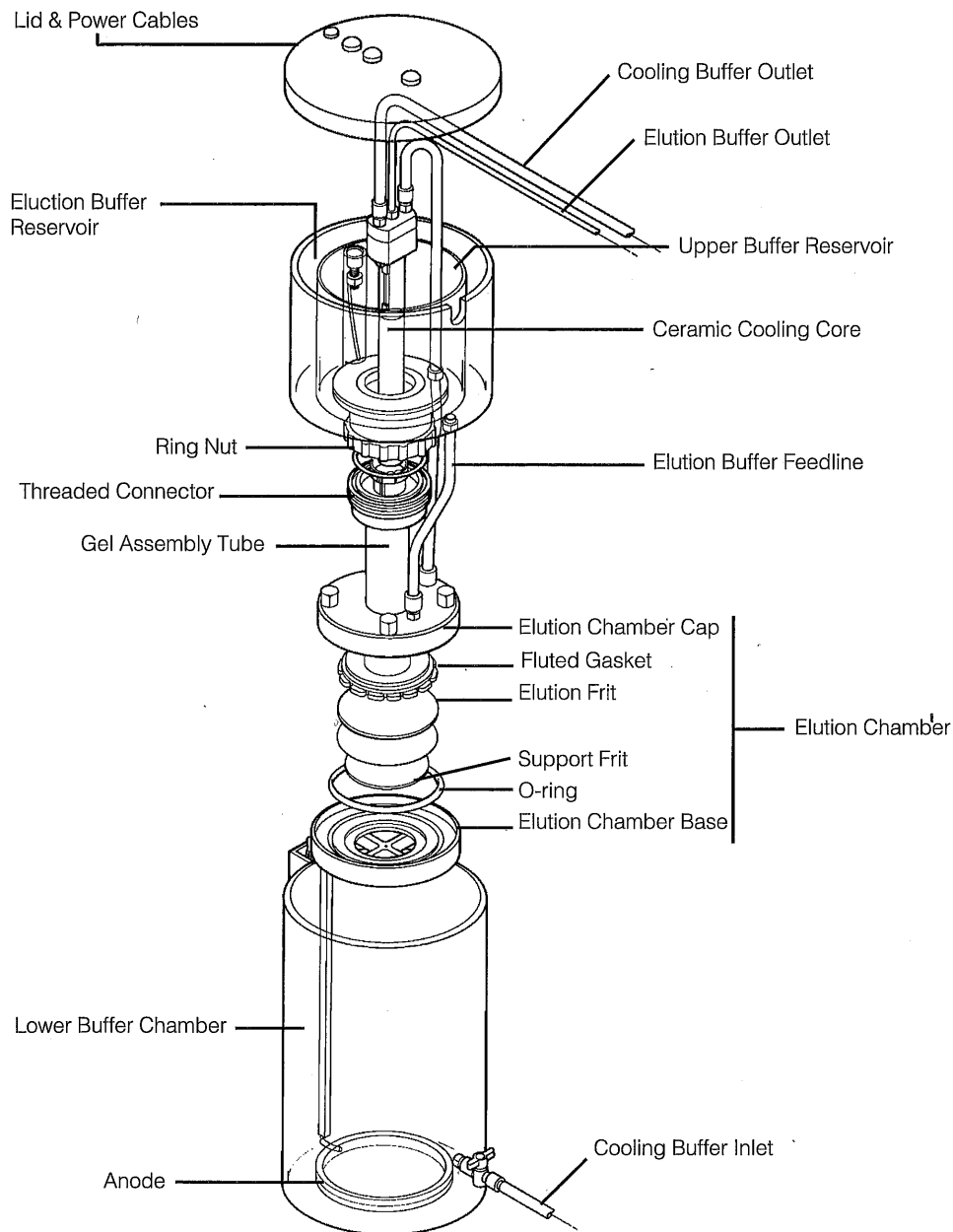
Description and Assembly of Major Components



Mini Prep Cell - Exploded View

Model 491 Prep Cell

Description of Major Components



Exploded View of the Model 491 Prep Cell

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

Website www.bio-rad.com U.S. (800) 4BIORAD Australia 02-9914-2800 Austria (1)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
 China (86-10) 2046622 Denmark 39 17 9947 Finland 90 804 2200 France (1) 43 90 46 90 Germany 089 31884-0 Hong Kong 7893300
 India 91-11-461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 The Netherlands 0313 18-540666 New Zealand 09-443 3099
 Singapore (65) 272-9877 Spain (91) 661 70 85 Sweden 46 (0) 8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800 181134