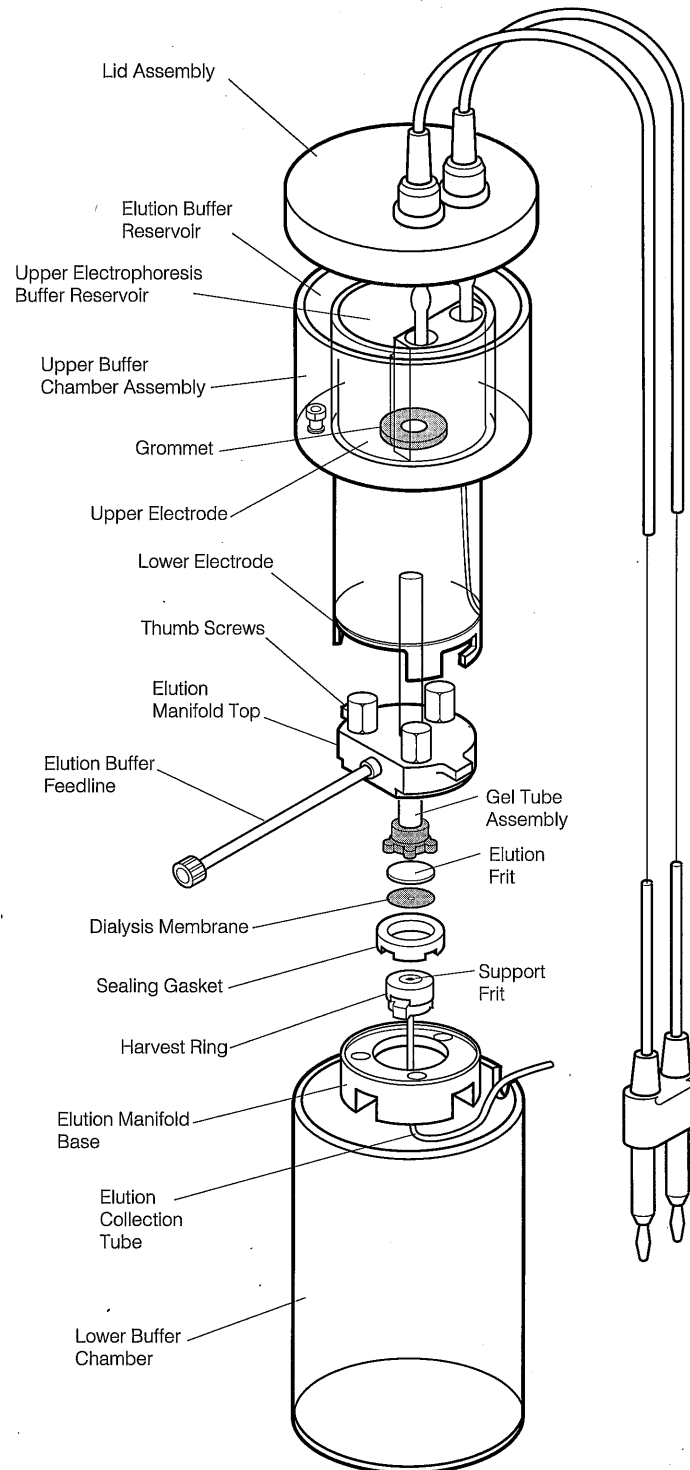




Mini Prep Cell

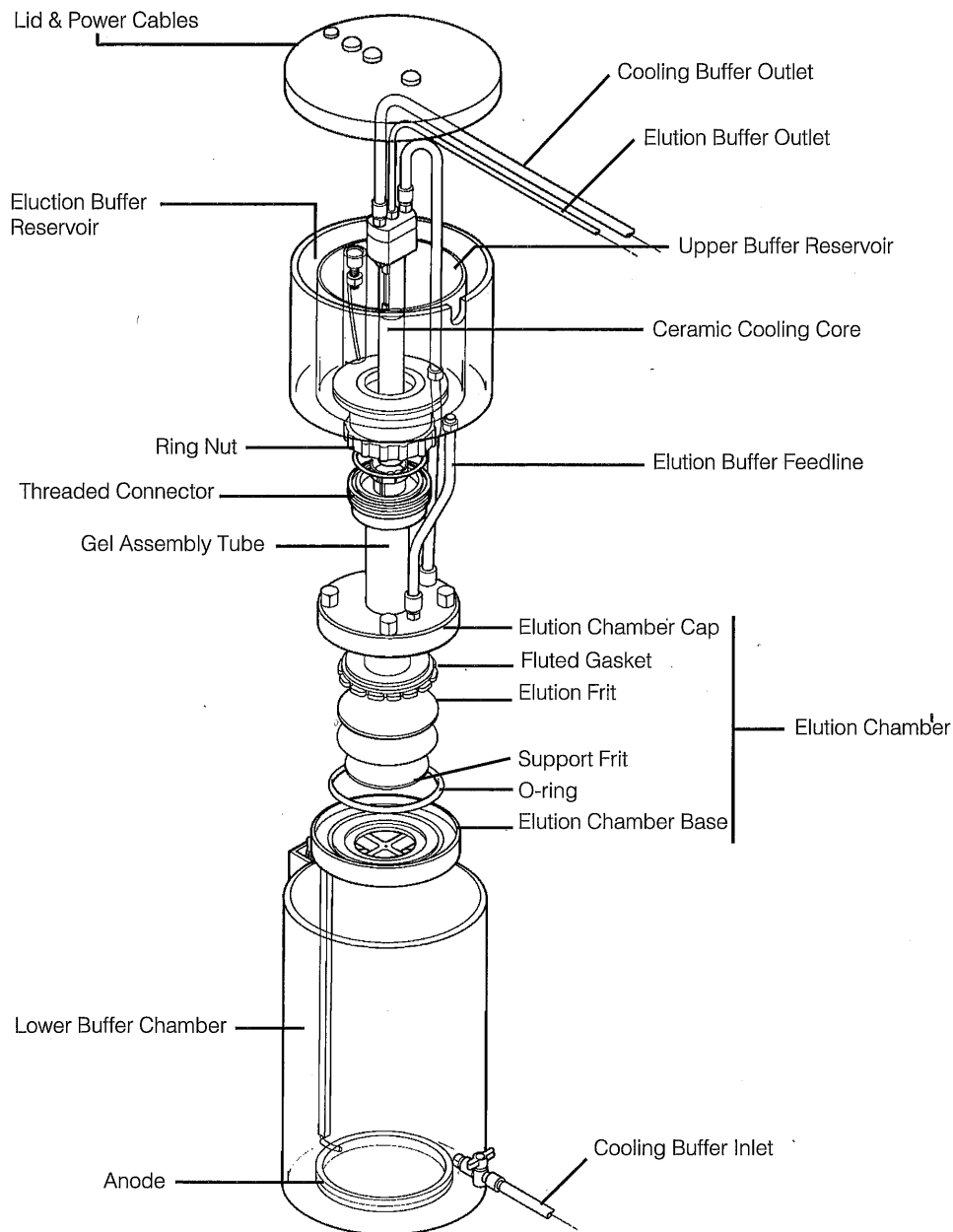
Description and Assembly of Major Components



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