



Mini Prep Cell and Model 491 Prep Cell Starter Kit Instructions

This is your Prep Cell starter kit. The kit has been provided to demonstrate how well the Prep Cell resolves two adjacent proteins and familiarize you with the assembly and operation of the Prep Cell. The sample in the starter kit consists of two prestained proteins, Soybean Trypsin Inhibitor (apparent MW 27,500 +/- 10%) and Carbonic Anhydrase (apparent MW 32,500 +/- 10%). Migration of these two blue proteins will demonstrate how straight these bands will migrate during the run. This run can be monitored. It is important to note that as a consequence of the staining process, the bandwidths of the prestained proteins are greater than those of unstained proteins.

Gel composition: 10% T/2.67% C - Resolving gel
4% T/2.67% C - Stacking gel

Gel length:	Mini Prep Cell	Model 491 Prep Cell
Resolving gel:	5.5 cm (2 ml)	5.5 cm (20 ml)
Stacking gel:	0.5 cm (200 µl)	0.5 cm (2 ml)
Gel tube size:	7 mm ID	28 mm ID
Sample load:	30 µl	300 µl (~150 µg of each protein)

All reagents necessary for three trial runs are included in this kit. It is advisable to perform at least one trial run to familiarize yourself with assembly and operation of the Prep Cell. Detailed diagrams and instructions are provided in the Prep Cell instruction manual. Please read the manual before your first run.

Reagent Preparation, Assembly, and Operation

Step 1: Prepare 30% T/2.67% C acrylamide monomer stock solution (100 ml)

0.8 g bis

29.2 g acrylamide

Dissolve in 70 ml distilled water. Add distilled water to a final volume of 100 ml. Filter and store at 4 °C (30 days maximum).

Step 2: Prepare 4x resolving gel buffer - 1.5M Tris/HCl, pH 8.8 (100 ml)

18.17 g Tris base

Dissolve in ~ 80 ml distilled water.

Adjust to pH 8.8 with 10 N HCl (DO NOT back titrate with base). Add distilled water to a final volume of 100 ml. Store at room temperature.

Step 3: Prepare 4x stacking gel buffer - 0.5M Tris/HCl, pH 6.8 (100 ml)

6.0 g Tris base

Dissolve in ~80 ml distilled water.

Adjust to pH 6.8 with 10 N HCl. (DO NOT back titrate with base).

Add distilled water to a final volume of 100 ml.

Store at room temperature.

Step 4: Prepare electrophoresis/elution buffer - Tris/Glycine/SDS (10 liters) Dilute 1 liter of 10x Tris/Glycine/SDS buffer with 9 liters of distilled water.

Step 5: Secure the gel tube assembly and cooling core to the casting stand and set up the cooling path needed during polymerization of the gel. See the instruction manual for details.

Step 6: Prepare resolving gel - 10% T/2.67% C, 0.375 M Tris/HCl, pH 8.8.

30% T/2.67% C stock solution 6.67 ml

1.5 M Tris/HCl, pH 8.8 5.0 ml

Distilled water 8.27 ml

10 % ammonium persulfate 50 µl*

TEMED 5 µl*

TOTAL MONOMER 20 ml

* Degas monomer solution prior to adding catalyst.

- Step 7:** Pour the degassed monomer solution with the added catalyst into the gel tube assembly. Carefully overlay with water or water saturated 2-butanol or tert-amyl alcohol. After ~ 2 hours replace overlay with 0.375 M Tris/HCl, pH 8.8 buffer and continue polymerization overnight. See the instruction manual for details.
- Step 8:** Prepare stacking gel - 4% T/2.67% C, 0.125 M Tris/HCl, pH 6.8
- | | |
|------------------------------|---------|
| 30% T/2.67% C stock solution | 1.33 ml |
| 0.5 M Tris/HCl, pH 6.8 | 2.5 ml |
| Distilled water | 6.1 ml |
| 10 % ammonium persulfate | 50 µl* |
| TEMED | 10 µl* |
| TOTAL MONOMER | 10 ml |
- * Degas monomer solution prior to adding catalyst.
- Step 9:** Carefully decant or aspirate the overlay buffer and add the stacking gel monomer solution. Carefully overlay with water, water saturated 2-butanol, or tert-amyl alcohol. Allow the gel to polymerize for 1–2 hours.
- Step 10:** Carefully decant or aspirate overlay. Assemble the elution chamber and the Prep Cell as described in the instruction manual.
- Step 11:** Add electrophoresis buffer to the lower and upper buffer chambers and to the elution buffer reservoir.
- Step 12:** Apply sample. Heat the sample to 40 °C for 1 minute to dissolve any solids which may have precipitated during storage. Using the sample application syringe, load the prestained proteins directly on the surface of the stacking gel, 300 µl for Model 491 Prep Cell and 30 µl for the Mini Prep Cell.

Recommended running conditions and expected results:

	Mini Prep Cell	Model 491 Prep Cell	
Power conditions:	4–5 mA constant, or 1 W constant	40 mA constant or 10–12 W constant	<ul style="list-style-type: none"> • The ion front and the free dye should elute in 2–2.5 hours. • Prestained Soybean Trypsin Inhibitor should elute after 2.5–3 hours, and prestained Carbonic Anhydrase after 3–3.5 hours. • Total electrophoresis time is approximately 4 hours.
Elution flow rate:	0.1 ml/min	1 ml/min	
Cooling flow rate:	N/A	80–100 ml/min	
Fraction collector:	0.25 ml/fraction	2.5 ml/fraction	

Product Information

Catalog

Number	Description
161-5104	Mini and Model 491 Prep Cell Starter Kit with Standard
161-5101	Mini and Model 491 Prep Cell Starter Kit without Standard
161-0323	Mini and Model 491 Prep Cell Standard



Bio-Rad
Laboratories

Life Science
Group

Bio-Rad Laboratories Main Office, 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-5800
 Also in: Reagents Park, Australia, Ph. 02-9914-2800, Fx. 02-9914-2888 Wien, Austria, Ph. (1) 877 89 01, Fx. (1) 876 56 29 Nazareth, Belgium, Ph. 09-385 55 11, Fx. 09-385 65 54
 Mississauga, Canada, Ph. (905) 712-2771, Fx. (905) 712-2990 Beijing, China, Ph. (01) 2046622, Fx. (01) 2051876 Copenhagen, Denmark, Ph. 39 17 9947, Fx. 39 27 1698
 Espoo, Finland, Ph. 90 804 2200, Fx. 90 804 1100 Ivry sur Seine Cedex, France, Ph. (1) 49 60 68 34, Fx. (1) 46 71 24 67 München, Germany, Ph. 089 31884-0, Fx. 089 31884-100
 New Delhi, India, Ph. 91-11-461-0103, Fx. 91-11-461-0765 Milano, Italy, Ph. 02-21609.1, Fx. 02-21609.399 Tokyo, Japan, Ph. 03-5811-6270 Fx. 03-5811-6272 Veendam,
 The Netherlands, Ph. 0318-540666, Fx. 0318-542216 Auckland, New Zealand, Ph. 09-443 3099, Fx. 09-443 3097 Kowloon, Hong Kong, Ph. 7893300, Fx. 7891257 Singapore,
 Ph. (65) 272-9877, Fx. (65) 273-4835 Solna, Sweden, Ph. 46 (0) 8 735 83 00, Fx 46 (0) 735 54 60 Madrid, Spain, Ph. (91) 661 70 85, Fx. (91) 661 96 98 Glattbrugg, Switzerland,
 Ph. 01/809 55 55, Fx. 01/809 55 00 Hemel Hempstead, United Kingdom, Ph. 0800 181134, Fx. 01442 259118