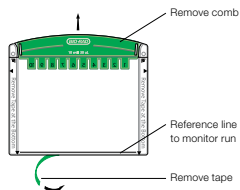


Instructions for Using Mini-PROTEAN Precast Gels

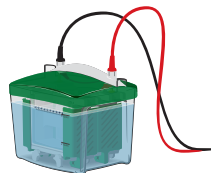
- 1. Remove Comb:** Position both thumbs on the ridges on the comb. Remove the comb by pushing upward in one smooth continuous motion.
- 2. Remove Tape:** Pull the green tape gently to remove it from the bottom of the cassette.
- 3. Assemble Mini-PROTEAN Tetra Cell:** Assemble the cassette into the running module of the Mini-PROTEAN Tetra system. Add running buffer to the inner and outer chambers. Use a syringe or a disposable transfer pipet to rinse the wells with running buffer.
- 4. Run Gel:** Prepare the samples and load into the wells. Run the gel until the dye front reaches the reference line. Refer to the instruction manual (Bulletin 1658100) for more information on running conditions. At the completion of the run, disconnect the cell and remove the cassette.
- 5. Open Cassette:** Align the arrow on the opening lever with the arrows marked on the cassette. Insert the lever between the cassette plates at all four locations and apply downward pressure to break the seal. Gently pull apart the two plates beginning from the top of the cassette.
- 6. Remove Gel:** Gently remove the gel from the cassette.

Please note: when running 1–2 gels in the Mini-PROTEAN Tetra cell, do not leave the companion module in the tank; do not run different gel types (chemistries) or percentages at the same time.

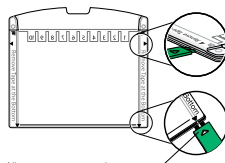
Preparing the Cassette



Running the Gel



Opening the Cassette



Align arrows on opening lever and cassette

Ordering Information

Description	 10-well 30 µl	 10-well 50 µl	 12-well 20 µl	 15-well 15 µl
Mini-PROTEAN Tris-Tricine Gels (Pack of 2)				
16.5% Resolving Gel	456-3063	456-3064	456-3065*	456-3066
10–20% Resolving Gel	456-3113	456-3114	456-3115*	456-3116*
Mini-PROTEAN TBE Gels (Pack of 2)				
5% TBE Gel	456-5013	456-5014*	456-5015*	456-5016
10% TBE Gel	456-5033	456-5034*	456-5035	456-5036
15% TBE Gel	456-5053*	456-5054	456-5055*	456-5056
4–20% TBE Gel	456-5093*	456-5094*	456-5095*	456-5096*
Mini-PROTEAN TBE-Urea Gels (Pack of 2)				
10% TBE-Urea Gel	456-6033*	—	—	456-6036*
15% TBE-Urea Gel	456-6053*	—	456-6055*	456-6056*

* Please allow up to 2 weeks for delivery.

Refer to the instruction manual (Bulletin 1658100) for more information on running conditions. Visit us at www.miniprotean.com for more information.

BIO-RAD