

Total Protein Staining

General protocols are described below for Mini-PROTEAN® gels. For more details, refer to the instruction manual for the stain you are using.

Bio-Safe Coomassie Stain

- 1 Wash gels three times for 5 min each in 200 ml diH₂O per gel.
- 2 Remove all water from staining container and add 50 ml of Bio-Safe Coomassie stain (or enough to completely cover gel). Agitate for 1 hr.
- 3 Rinse in 200 ml diH₂O for ≥30 min. Stained gels can be stored in water.

Oriole Fluorescent Gel Stain

- 1 If using the 5 L configuration, prepare Oriole stain solution by adding 400 ml of methanol to the 1 L bottle of diluents. Then add 10 ml of Oriole fluorescent gel stain concentrate and mix well by shaking.
- 2 Place gel in a staining tray with 50 ml of Oriole fluorescent gel stain. Cover the tray, place on a rocker, and agitate gently for ~1.5 hr.
- 3 Transfer the gel to diH₂O prior to imaging. Destaining is not necessary.

Flamingo Fluorescent Gel Stain

- 1 Place gel in a staining tray with 100 ml of fixing solution (40% ethanol, 10% acetic acid). Cover the tray, place on a rocker, and agitate gently for at least 2 hr.
- 2 Pour off the fix solution and add 50 ml of 1x stain solution (dilute 1 part Flamingo fluorescent gel stain with 9 parts diH₂O). Cover the tray, place on a rocker or shaker and agitate gently. Stain for at least 3 hr.
- 3 (Optional) Carefully pour off the stain solution and replace with an equal volume of 0.1% (w/v) Tween 20. Cover the tray, place on a rocker or shaker and agitate gently for 10 min.
- 4 Rinse gel with diH₂O prior to imaging.

TIPS

- Always wear gloves during the staining process. Try to avoid touching the gels with your fingers. Wet gloves with water or buffer before handling the gel to keep the gel from sticking and tearing
- Use clean and dust-free containers for gel staining. If possible, place a lid on the container to avoid contamination of the staining solution
- Gently agitate the container on a horizontal shaker, making sure the gel is completely covered with stain solution all the time
- Use pure chemicals and highly purified diH₂O (conductivity <2 μS)
- Fluorescent dyes like Flamingo and Oriole fluorescent gel stains have a higher dynamic range than Coomassie or silver staining techniques, making them suitable for quantitative protein analysis
- Gels stained with fluorescent dyes can be counterstained with colloidal Coomassie for further reference. In fact, doing so enhances sensitivity of the colloidal Coomassie stain
- For long term-storage, shrink-wrap the stained gels in a 10% glycerol solution (storage at 4°C) or dry them with a GelAir™ drying system within 60 min

Silver Staining (Bio-Rad Silver Stain)

Step	Reagent	Volume	Duration		
			<0.5 mm Gel	0.5–1.0 mm Gel	>1.0 mm Gel
1	Fixative 40% methanol, 10% acetic acid	400 ml	30 min	30 min	60 min
2	Fixative 10% ethanol/5% acetic acid	400 ml	15 min	15 min	130 min
3		400 ml	15 min	15 min	130 min
4	Oxidizer	200 ml	3 min	5 min	10 min
5	diH ₂ O	400 ml	2 min	5 min	10 min
6		400 ml	2 min	5 min	10 min
7	(Repeat washes 5–7 until all the yellow color is gone from the gel)	400 ml	2 min	5 min	10 min
8	Silver reagent	200 ml	15 min	20 min	30 min
9	diH ₂ O	400 ml	—	1 min	2 min
10	Developer	200 ml	~30 sec. Develop until solution turns yellow or until brown precipitate appears.		
11		200 ml	~5 min	~5 min	~5 min
12		200 ml	—	~5 min	~5 min
13	Stop 5% acetic acid	400 ml	~5 min	~5 min	~5 min

This is an excerpt from Bio-Rad's comprehensive Electrophoresis Guide (Bulletin 6040).



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000