

Western Blotting Protocol: Mini-PROTEAN TGX Stain-Free Gel Run on a Tank Transfer System Using 1x Tris Buffered Saline (TBS) with 1% Casein and Fluorescence Detection Protocol

Introduction

This easy protocol includes acceptable ranges for each applicable step. This protocol can be printed and used at the bench as a checklist for tracking the actual conditions used.

Electrophoresis

✓ Step	Instructions	Actual Conditions	Low End of Range	Ideal Conditions	High End of Range
1	Add 4x or 2x Laemmli buffer to the sample for a final concentration of 1x. Prepare 50% more than you intend to load.				
2	Heat at 95°C for 5 min.		70 2	95°C 5 min	100 10
3	Centrifuge at 10,000 x g for 5 min.		5,000 2	10,000 x g 5 min	16,000 30
4	Insert the gel and fill the gel apparatus to fill line (550 ml for one to two gels and 800 ml for three to four gels) with running buffer.				
5	Load 10 µl protein ladder/molecular weight standard.		1	10 µl	25
6	Load the gel with desired samples.				
7	Run the gel at 300 V until the dye reaches the bottom of the gel.		100	300 V	350
8	Remove the gel from the cassette and place on the imaging tray.				
9	Activate the Stain-Free dye and acquire a gel image.				

Transfer

✓ Step	Instructions	Actual Conditions	Low End of Range	Ideal Conditions	High End of Range
1	Equilibrate the transfer buffer to the appropriate temperature.		4	25°C	30
2	Equilibrate the membrane, filter paper, and fiber pads in transfer buffer for 20 min.				
3	Prepare the gel sandwich and place the cassette into the module.				
4	Add cooling unit, stir bar, and transfer buffer.				
5	Run the transfer at 100 V for 1 hr.		5 0.5	100 V 1 hr	200 16
6	Rinse the membrane with 1x TBS and place on the blot tray.				
7	Acquire a blot image.				

Blocking

✓	Step	Instructions	Actual Conditions	Low End of Range	Ideal Conditions	High End of Range
	1	Immerse the membrane in 10 ml of 1x TBS with 1% casein.				
	2	Place on a rocker and block at room temperature (RT) for 1 hr.		0.5	1 hr	3

Primary Immunodetection

✓	Step	Instructions	Actual Conditions	Low End of Range	Ideal Conditions	High End of Range
	1	Dilute the antibody/antibodies in 10 ml of 1x TBS with 1% casein. Dilution ratio: 1: _____				
	2	Place on a rocker and incubate at RT for 1 hr.		4°C 0.5	RT 1 hr	30°C 12
	3	Wash with 10 ml of 1x TBS with 0.05% Tween 20 for 5 min with agitation.		0.01 3	1x TBS with 0.05% Tween 20 5 min	0.1 10
	4	Repeat step 3 for a total of five washes.		3	5 cycles	7

Secondary Immunodetection

✓	Step	Instructions	Actual Conditions	Low End of Range	Ideal Conditions	High End of Range
	1	Dilute the antibody/antibodies in 10 ml of 1x TBS with 0.05% Tween 20. Dilution ratio: 1: _____				
	2	Place on a rocker and incubate at RT for 1 hr.		4°C 0.5	RT 1 hr	30°C 12
	3	Wash with 10 ml of 1x TBS with 0.05% Tween 20 for 5 min with agitation.		0.01 3	1x TBS with 0.05% Tween 20 5 min	0.1 10
	4	Repeat step 3 for a total of six washes.		4	6 cycles	8

Detection

✓	Step	Instructions	Actual Conditions	Low End of Range	Ideal Conditions	High End of Range
	1	Acquire a series of images until saturation of the target band(s) is observed.				
	2	Acquire a final image with no saturated bands. Filename: _____				

Signed _____ Date _____

Countersigned _____ Date _____

Go to bio-rad.com/WesternResources for more information, tips, tricks, and troubleshooting.

BIO-RAD and MINI-PROTEAN are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

TGX Stain-Free Precast Gels are covered by U.S. Patent Numbers 7,569,130 and 8,007,646.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 6511 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

