

Western Blotting Protocol: Mini-PROTEAN TGX Stain-Free Gel Run on a Tank Transfer System Using EveryBlot Blocking Buffer and Fluorescence Detection

Protocol

Western Blotting

Bulletin 7440

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 Tris buffered saline (TBS) and place on the blot tray. | | | | |
| 7 | Acquire a blot image. | | | | |

BIO-RAD

Blocking

| ✓ Step | Instructions | Actual Conditions | Low End of Range | Ideal Conditions | High End of Range |
|--------|---|-------------------|------------------|------------------|-------------------|
| 1 | Immerse the membrane in 10 ml of EveryBlot Blocking Buffer. | | | | |
| 2 | Place on a rocker and block at room temperature (RT) for 5 min. | | 3 | 5 min | 30 |

Primary Immunodetection

| ✓ Step | Instructions | Actual Conditions | Low End of Range | Ideal Conditions | High End of Range |
|--------|---|-------------------|------------------|----------------------------------|-------------------|
| 1 | Dilute the antibody in 10 ml of EveryBlot Blocking Buffer. 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 TBS with 0.05% Tween 20 for 5 min with agitation. | | 0.01 3 | 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 in 10 ml of EveryBlot Blocking Buffer. 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 TBS with 0.05% Tween 20 for 5 min with agitation. | | 0.01 3 | 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 |
|--------|--|
| 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 2 651 8311
United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

