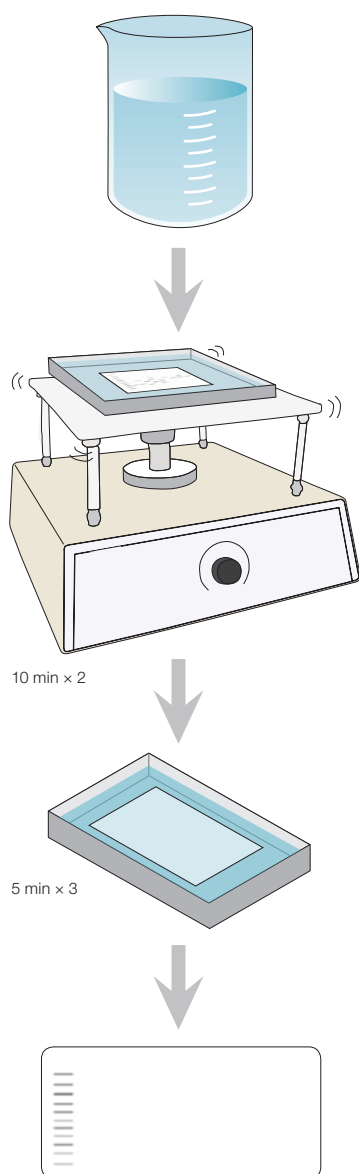


# Blot Stripping and Reprobing

Protocol

Bulletin 6218

## Based on Legocki and Verma 1981



1

Prepare acidic glycine stripping buffer (0.1 M glycine, 20 mM magnesium acetate, 50 mM KCl, pH 2.2; for recipe).

2

Add enough acidic glycine stripping buffer to completely cover the developed membrane and incubate at room temperature for 10 min with gentle agitation.

3

Repeat step 2 with fresh acidic glycine stripping buffer.

4

Wash the blot three times in TTBS for 5 min each with gentle agitation.

5

Test for complete removal of primary antibody by reprobing with only the secondary antibody and redeveloping. No signal should be detectable. Re-block the membrane and proceed to the next detection protocol.

### TIPS

This protocol (based on Legocki and Verma 1981) uses low pH to gently remove antibody from the membrane. The protocol removes little of the sample proteins but may not remove all antibodies with high affinities for their targets

**BIO-RAD**

This is an excerpt from Bio-Rad's comprehensive Protein Blotting Guide (Bulletin 2895).



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

---

**Web site** [www.bio-rad.com](http://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

---