



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Western ReProbe™ PLUS

For Stripping High Affinity Antibodies
from Western Blots

(Cat. # 786-307, 786-308, 786-309)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

Based on our popular Western ReProbe™, the modified formulation allows for the removal of stubborn, high affinity antibodies from membrane bound proteins without destroying the antigenic binding affinity. The membrane bound protein is retained on the membrane and the matching antibodies are washed away. Once the antigen-antibody bonds are broken, the membrane bound protein is free to accept new antibodies. Each 100ml of Western ReProbe™ PLUS is sufficient for 25-30 standard size (7.5cm x 8.5cm) western blots.

Western ReProbe™ PLUS is not recommended for stripping color producing Western blots that use substrates such as TMB, chloronaphthol and DAB.

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-307	Western ReProbe™ PLUS	500ml
786-308	Western ReProbe™ PLUS	1L
786-309	Western ReProbe™ PLUS	1gal

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store at room temperature. Use sterile conditions for removing solution from the bottle. When used properly, it is stable for 1 year.

ADDITIONAL ITEMS REQUIRED

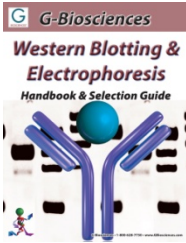
Reagents for blocking the membrane and/or immunodetection reagents

PROTOCOL

1. Following probing and detection of antibodies keep the membrane in an appropriate wash buffer (i.e. PBS + 0.1% Tween® (PBST) or TBS + 0.1% Tween® (TBST)).
2. Incubate the membrane in undiluted Western ReProbe™ PLUS buffer for 30 minutes at room temperature while shaking. Increasing the incubation time to 45 minutes would improve membrane stripping.
3. Wash the membrane 3 times with TBST or PBST. Use a large volume of washing buffer in a tray for washing. Each wash cycle should consist of agitating the membrane for 5 minutes in the washing buffer.
4. Block the membrane and perform immunodetection using normal protocols.
NOTE: Western ReProbe™ PLUS can be used for stripping and reprobing membranes 2-3 times. Some loss of signal may be noticed after the second cycle of stripping and re-probing.

RELATED PRODUCTS

Download our Western Blotting Handbook.



<http://info.gbiosciences.com/complete-western-blot-handbook--selection-guide/>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 8/3/2012 CMH



www.GBiosciences.com