



459PR-01

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Western ReProbe™

## For Stripping and Reprobing Western Blots

(Cat. # 786-119, 786-119S, 786-305, 786-306, 786-818)



**think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)**

## INTRODUCTION

Western ReProbe™ breaks antigen-antibody 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 proteins. Each 100ml of 1X Western ReProbe™ is sufficient for 25-30 standard size (7.5cm x 8.5cm) western blots. Western ReProbe™ is not recommended for stripping color producing Western blots that use substrates such as TMB, chloronaphthol and DAB.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-119	Western ReProbe™ [5X]	100ml
786-119S	Western ReProbe™ [5X]	10ml
786-305	Western ReProbe™ [5X]	500ml
786-306	Western ReProbe™ [5X]	1L
786-818	Western ReProbe™ [5X]	1 Gallon

## STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store at 4°C. 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

## PREPARATION BEFORE USE

1. Prepare the required volume (15-20ml/ blot) of 1X Western ReProbe™ by diluting the supplied 5X Western ReProbe™ in DI water (e.g. take 5ml of 5X Western ReProbe™ and add to 20ml DI water to make 25ml of 1X Western ReProbe™).

## 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 1X Western ReProbe™ 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.

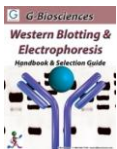
***NOTE:*** Western ReProbe™ can be used for stripping and reprobing membranes 2-3 times. Some loss of signal may be noticed after the second cycle of stripping and re-probing.

## CITATIONS

1. Ebrahem, Q. et al (2011) Invest. Ophthalmol. Vis. Sci. 52:6117
2. George, J. et al (2010) Neuro Oncology. doi:10.1093/neuonc/noq079
3. Qi, J. et al (2009) J. Biol. Chem. 284: 19927 - 19936
4. George, J. et al (2009) Clin. Cancer Res. 15: 7186 - 7195
5. Di, X. et al (2008) Hum. Reprod. 23: 1873-83
6. Roy, S. and Tenniswood, M. (2007) J. Biol. Chem. 282: 4765
7. George, J. et al (2007) Clin Cancer Res. 13(12): 3507-17
8. Joshi, M. et al (2006) Amer J. Respir. 10.1164
9. Brown, K. et al (2006) Am. J. Physiol. Lung 290: 259
10. Bu, S. et al (2006) Reproduction 131:1099
11. Lam, A.M.I and Frick, D.N. (2006) J. Virol. 80: 404
12. Roy, S. et al (2005) Cell Death Differ. 12: 482
13. Li, Q. et al (2004) Reproduction 128:555
14. Bakke, L.J. et al (2004) Biol. Reprod. 71:605
15. Rubin, M. et al (2004) Cancer Res. 64: 3814
16. Small, G. et al (2004) Mol. Pharmacol. 66: 1478
17. Small, G. et al (2003) J. Pharmacol. Exp. Ther. 307: 861
18. Kobori, H. et al (2003) Hypertension 41: 592
19. Kobori, H. et al (2003) Hypertension 41: 42
20. Kuefer, S. et al (2002) Am. J. Path. 161: 841
21. Muthumani, K. et al (2002) JBC 277: 37820
22. Dash, A. et al (2002) Am. J. Path. 161: 1743
23. Orlawski, R. et al (2002) J. Biol. Chem. 277: 27864
24. Schraufstatter, I. et al (2002) J. Immunol. 169: 2102
25. Schraufstatter, I. et al (2002) J. Immunol. 171: 6714
26. Zhang, Y. et al (2001) Mol. Endocrinol. 15: 1891

## 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](http://www.GBiosciences.com) or contact us.



[www.GBiosciences.com](http://www.GBiosciences.com)