

# Restore™ Plus Western Blot Stripping Buffer

46428 46430 46431

1959.4

Number	Description
46428	Restore Plus Western Blot Stripping Buffer, 30mL
46430	Restore Plus Western Blot Stripping Buffer, 500mL
46431	Restore Plus Western Blot Stripping Buffer, 5L

**Storage:** Upon receipt store at room temperature. Product shipped at ambient temperature.

## Introduction

Thermo Scientific Restore Plus Western Blot Stripping Buffer provides a robust but gentle method for removing primary and secondary antibodies from Western blots that were detected with chemiluminescent substrates. Restore Plus Western Blot Stripping Buffer allows reprobing saving time and costs when samples are in limited quantities, when the same sample requires analysis by different antibodies and when optimization is required. Traditional stripping methods either adversely alter the proteins on the membrane or use conditions that are effective for only low-affinity antibody-antigen interactions.<sup>1,2</sup> Restore Plus Western Blot Stripping Buffer can be used at room temperature to quickly and effectively strip most high-affinity antigen-antibody interactions.

## Additional Materials Required

- Western blot, previously blocked, probed and detected with a chemiluminescent substrate
- Wash Buffer such as Tris-buffered saline (TBS, Product No. 28376) or phosphate-buffered saline (PBS, Product No. 28372) containing 0.05% Tween®-20 (Product No. 28320)
- Primary and secondary antibodies
- Film or CCD camera for detecting the chemiluminescent signal

## Protocol for Stripping an Immunoblot

### Notes:

- When performing multiple strips and reprobing for different antigens, probe for the low-abundant proteins first.
  - Restore Plus Western Blot Stripping Buffer will not dissociate interactions between a biotinylated target protein and avidin-conjugated probes.
  - Stripping and reprobing fluorescent Western blots is not recommended because results are typically inconsistent.
1. Wash blots in Wash Buffer to remove the chemiluminescent substrate. Blots may be stored in PBS or TBS at 4°C until the stripping procedure can be performed.
  2. Place the blot in Restore Plus Western Blot Stripping Buffer and incubate for 5-15 minutes at room temperature. Use a sufficient volume to ensure that the blot is completely wetted (i.e. approximately 20mL for an 8 × 10cm blot).  
**Note:** Optimization of both incubation time and temperature is essential for best results. In general, high-affinity antibodies require at least 15 minutes and may require incubation at 37°C.
  3. Remove the blot from the Restore Plus Western Blot Stripping Buffer and wash in Wash Buffer.
  4. Block membrane for 30-60 minutes at room temperature or overnight at 2-8°C.

5. Test for removal of the antibodies as follows:
  - To test for complete removal of the enzyme conjugate (e.g., secondary antibody), incubate the membrane with new chemiluminescent substrate working solution and expose blot to film. If no signal is detected using a minimum 5 minute exposure, the enzyme conjugate has been successfully removed from the antigen or primary antibody.
  - To test for complete removal of the primary antibody, incubate the membrane with the enzyme conjugate and then wash in Wash Buffer. Incubate in new chemiluminescent substrate working solution and expose to film. If no signal is detected with a 5 minute exposure, the primary antibody has been successfully removed from the antigen.
6. If signal is detected with either test in Step 5, return to Step 1, stripping for an additional 5-15 minutes. Some antigen/antibody systems require increased temperature and/or longer incubation times. Optimize stripping time and temperature to ensure complete removal of antibodies while minimizing damage to the antigen.
7. After determining that the membrane is properly stripped, the second immunoprobng experiment may be performed.
 

**Note:** Blot may be stripped and reprobng several times but might require longer exposure times or a more sensitive chemiluminescent substrate. Subsequent reprobngs might result in decreased signal if the antigen is labile in Restore Plus Western Blot Stripping Buffer. Analysis of the individual system is required.

## Troubleshooting

Problem	Possible Cause	Solution
Background after stripping and subsequent detection	Not sufficiently blocked after stripping	Optimize blocking conditions
	Antibody concentrations are too high	Strip again and reprobe using more dilute antibody concentrations
Loss of signal or no signal after stripping and reprobng	Antigen is not present or in low abundance	Prepare a new blot and probe for low-abundant antigens first
	Protein or antibody concentrations were too low	Increase antibody concentrations
		Load more protein in the gel
Signal obtained after stripping	Extremely high-affinity antigen-antibody interaction	Increase incubation time and temperature

## Related Thermo Scientific Products

- 46640** SuperSignal Western Blot Enhancer, 500mL kit, sufficient for 25 mini blots or 2000cm<sup>2</sup>
- 46641** SuperSignal Western Blot Enhancer, 50mL kit, sufficient for 2 mini blots or 160cm<sup>2</sup>
- 32106** Pierce<sup>®</sup> ECL Western Blotting Substrate, 500mL
- 34080** SuperSignal<sup>®</sup> West Pico Chemiluminescent Substrate, 500mL
- 34075** SuperSignal West Dura Chemiluminescent Substrate, 100mL
- 34095** SuperSignal West Femto Chemiluminescent Substrate, 100mL

## General References

1. Kaufmann, S.H., *et al.* (1987). The erasable Western blot. *Anal Biochem* **161**: 89-95.
2. Kaufmann, S.H. and Kellner, U. (1998). Erasure of Western blots after autoradiographic or chemiluminescent detection. In *Immunochemical Protocols*. Ed. Pound, J.D. Humanna Press, Totowa, NJ., p. 223-35.

Tween is a trademark of Croda International Plc.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

**No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).**

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.