

Restore Western Blot Stripping Buffer

21059 21062 21063

0887.6

Number	Description
21062	Restore Western Blot Stripping Buffer, 30mL
21059	Restore Western Blot Stripping Buffer, 500mL
21063	Restore Western Blot Stripping Buffer, 5L

Storage: Upon receipt store at room temperature or 4°C. Product shipped at ambient temperature.

Introduction

Nitrocellulose and PVDF membranes probed by Western blotting procedures and detected by chemiluminescent or other non-precipitating substrates can be stripped and re-probed using Thermo Scientific™ Restore™ Western Blot Stripping Buffer. Western blotting is widely used to detect and compare proteins in complex mixtures, and chemiluminescence has largely replaced chromogenic substrates as the most convenient and sensitive method of detection. One advantage of chemiluminescence is the ability to strip and reprobe the protein mixture on the membrane. Traditional stripping methods use conditions that are effective for only low-affinity antibody-antigen interactions or are so harsh that they tend to adversely alter the antigen for subsequent immunoprobings.^{1,2} Restore Western Blot Stripping Buffer is a robust but gentle formulation for stripping primary and secondary antibodies from blots to enable several reprobings on the same membrane. Restore Western Blot Stripping Buffer is ideal for use with Thermo Scientific™ SuperSignal™ West Chemiluminescent Substrates.

Additional Materials Required

- Western blot, previously blocked, probed and detected with chemiluminescent (i.e., nonprecipitating) substrate
- Wash buffer such as Tris-buffered saline (TBS) or phosphate-buffered saline (PBS) with 0.05% Tween™ 20
- Primary and secondary antibodies for both first and second Western blotting experiments
- Film, the MYECL™ Imager or other imaging instrument for detecting the chemiluminescent signal

Procedure for Stripping an Immunoblot

Notes:

- Blots may be stored in PBS or TBS at 4°C until the stripping procedure can be performed.
- Restore 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. Warm the bottle of Restore Western Blot Stripping Buffer to room temperature.
2. Place the blot in Restore Western Blot Stripping Buffer and incubate for 5 to 15 minutes at 37°C. Use a sufficient volume to ensure that the blot is completely wetted (i.e., ~20mL required for an 8 × 10cm blot).

Note: Optimization of both incubation time and temperature is essential for best results. For some antibodies, room temperature incubation is sufficient. However, high-affinity antibodies or saturated blots (excess secondary antibody) may require incubation for an additional 5 to 10 minutes at 37°C.

3. Remove the blot from the Restore Western Blot Stripping Buffer and wash in TBS or PBS.

4. Test for the removal of the immunodetection reagents as follows:
 - Test for complete removal of the HRP label (e.g., secondary antibody): Incubate the membrane with new SuperSignal West Working Solution and expose to film. If no signal is detected using a 5-minute exposure, the HRP conjugate has been successfully removed from the antigen or primary antibody.
 - Test for complete removal of the primary antibody: Incubate the membrane with the HRP-labeled secondary antibody, followed by a wash in wash buffer. Incubate in new SuperSignal West 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.
5. If signal is detected with either test in step 4, return to step 2, stripping for an additional 5-15 minutes. Some antigen/antibody systems require increased temperature and/or longer incubation times to strip them fully. Optimize stripping time and temperature to ensure complete removal of antibodies while preventing damage to the antigen.
6. After determining that the membrane is properly stripped, the second immunoprobng experiment may be performed.

Notes:

- Blot can be stripped and reprobed several times but might require longer exposure times or a more sensitive chemiluminescent substrate. Subsequent reprobings might result in decreased signal if the antigen is labile in Restore Western Blot Stripping Buffer. Analysis of the individual system is required.
- Reblocking the membrane is recommended after stripping.

Related Thermo Scientific Products

62236	MYECL Imager
62288	Pierce™ G2 Fast Blotter
46640	SuperSignal Western Blot Enhancer, 500mL kit, sufficient for 25 mini blots or 2000cm²
34080	SuperSignal West Pico Chemiluminescent Substrate, 500mL
34075	SuperSignal West Dura Extended Duration Substrate, 100mL
34095	SuperSignal West Femto Maximum Sensitivity Substrate, 100mL
28374	BupH™ Modified Dulbecco's PBS Packs, 40 packs, each pack yields 500mL of buffer
28376	BupH Tris Buffered Saline Packs, 40 packs, each pack yields 500mL of buffer

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. Humana Press, Totowa, NJ., p. 223-35.

Products are 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"). 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 Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

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 REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR 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.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

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

© 2014 Thermo Fisher Scientific Inc. All rights reserved. Tween is a trademark of Croda International Plc. Unless otherwise indicated, all other trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.