


Restore™ Western Blot Stripping Buffer

Doc. Part No. 2160887 Pub. No. MAN0011403 Rev. B.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

Thermo Scientific™ Restore™ 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™ Western Blot Stripping Buffer allows for 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. Restore™ Western Blot Stripping Buffer can be used with both nitrocellulose and PVDF-probed membranes.

Contents and storage

Product	Cat. No.	Amount	Storage
Restore™ Western Blot Stripping Buffer	21062	30 mL	Room temperature or 4°C
	21059	500 mL	
	21063	5 L	

Procedural guidelines

- Restore™ Western Blot Stripping Buffer is supplied in ready-to-use format. It is not recommended to dilute the buffer.
- 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 fluorescent western blots is not recommended with Restore™ Western Blot Stripping Buffer as results can be inconsistent. It is recommended to use Restore™ Fluorescent Western Blot Stripping Buffer for fluorescent applications.
- When performing multiple strippings and reprobings for different antigens, it is recommended to probe for low-abundant proteins first.
- Reblocking the membrane is recommended after stripping.

Strip western blots and test effectiveness of procedure

1. Bring Restore™ Western Blot Stripping Buffer to room temperature.
2. Wash blot once for 5 minutes in wash buffer (PBS or TBS) to remove the chemiluminescent substrate.
3. Add sufficient Restore™ Western Blot Stripping Buffer to cover the blot and incubate for 5 to 15 minutes at 37 °C with gentle shaking.

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.

4. Remove the blot from the Restore™ Western Blot Stripping Buffer and wash 2 times for 5 minutes each in TBS or PBS with gentle shaking.

5. Test the effectiveness of stripping as follows:

To test for complete removal of the alkaline phosphatase (AP) or horseradish peroxidase (HRP) label (e.g., secondary antibody), incubate the membrane with new chemiluminescent substrate. If no signal is detected using a 5-minute exposure, the AP or HRP 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 AP or HRP-labeled secondary antibody, followed by 3, 5-minute washes in wash buffer. Incubate the membrane with new chemiluminescent substrate. 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, incubate blot in Restore™ Western Blot Stripping Buffer for an additional 5 to 15 minutes and repeat step 3 and step 4.

Note: 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.

7. After determining that the membrane is properly stripped, the second immunoprobng experiment may be performed.

8. Block the membrane in blocking buffer before continuing with the second immunoprobng experiment.

Related products

Table 1 Stripping buffers for western blotting

Product	Select when ...	Available formats
Restore™ Western Blot Stripping Buffer	Primary antibody is susceptible to stripping buffers	30 mL (Cat. No. 21062) 500 mL (Cat. No. 21059) 5 L (Cat. No. 21063)
Restore™ PLUS Western Blot Stripping Buffer	Removing high affinity antibodies	30 mL (Cat. No. 46428) 500 mL (Cat. No. 46430)
Restore™ Fluorescent Western Blot Stripping Buffer	Removing fluorescent-labeled antibodies	20 mL (Cat. No. 62299) 100 mL (Cat. No.62300)

Table 2 Additional products

Products	Learn more
ECL Kits	thermofisher.com/chemisubstrates
Western blot blocking buffers	thermofisher.com/blockingbuffers
Western blot transfer equipment and supplies	thermofisher.com/transfer
Western blot imaging and analysis	thermofisher.com/westernimaging

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0011403

Revision	Date	Description
B.0	13 December 2021	Updated content and move to CCMS
A.0	17 October 2015	New document

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

