

# SuperSignal<sup>®</sup> Western Blot Enhancer

46640 46641

2265.1

<b>Number</b>	<b>Description</b>
46640	<b>SuperSignal Western Blot Enhancer</b> , sufficient material for 25 mini blots or 2000cm <sup>2</sup> of membrane <b>Kit Contents:</b> <b>Antigen Pretreatment Solution</b> , 250mL <b>Primary Antibody Diluent</b> , 250mL
46641	<b>SuperSignal Western Blot Enhancer</b> , sufficient material for two mini blots or 160cm <sup>2</sup> of membrane <b>Kit Contents:</b> <b>Antigen Pretreatment Solution</b> , 25mL <b>Primary Antibody Diluent</b> , 25mL

**Storage:** Upon receipt store at 4°C. Product is shipped at ambient temperature.

## Introduction

The Thermo Scientific SuperSignal Western Blot Enhancer increases both signal intensity and sensitivity 3- to 10-fold compared to detection performed using conventional Western blotting. When a protein or antigen is difficult to detect because of low abundance or poor immunoreactivity, the SuperSignal Western Blot Enhancer can significantly reduce background and enhance detection of low abundant and weakly immunoreactive antigens. This versatile kit works with PVDF and nitrocellulose membranes and is compatible with fluorescence, chromogenic and chemiluminescent detection.

## Important Product Information

- Apply the Antigen Pretreatment Solution to the membrane after protein transfer and before applying blocking buffer.
- Use the Primary Antibody Diluent to dilute the primary antibody.
- Use any blocking solution to prepare appropriate dilution of secondary antibody conjugate.

## Additional Materials Required

- Ultrapure water
- Nitrocellulose or PVDF membrane with transferred proteins
- Appropriate primary antibody and secondary antibody conjugate
- Wash Buffer: phosphate- or Tris-buffered saline containing 0.05% Tween<sup>®</sup>-20 Detergent
- Blocking Buffer (e.g., Thermo Scientific SuperBlock Blocking Buffer, Product No. 37515 or 37535)
- Shaker for membrane incubations

## Procedure for Western Blot Enhancer Treatment

**Important Note:** For all steps, use sufficient volumes to completely immerse the membrane. Enhancer treatment must be completed before performing membrane blocking.

1. Perform electrophoresis and transfer protein onto a nitrocellulose or PVDF membrane.
2. Wash the membrane with ultrapure water for two minutes with shaking.
3. Discard the water and rinse the membrane twice with ultrapure water.
4. Add sufficient volume of Antigen Pretreatment Solution to immerse the membrane. Incubate at room temperature for 10 minutes with shaking.
5. Discard the solution from the tray. Rinse the membrane five times with ultrapure water.
6. Incubate the membrane with Blocking Buffer.
7. Rinse the membrane with Wash Buffer three times. Wash the membrane with Wash Buffer for 5 minutes with shaking.
8. Dilute the primary antibody in Primary Antibody Diluent. Use an antibody concentration appropriate for the specific detection method. See Table 1 for concentration ranges for chemiluminescent substrates.
9. Incubate the membrane in the diluted primary antibody for 1 hour at room temperature with shaking.
10. Wash the membrane with Wash Buffer four times for 5 minutes each.
11. Dilute the secondary antibody conjugate in Blocking Buffer. Use an antibody concentration appropriate for the specific detection method. See Table 1 for concentration ranges for chemiluminescent substrates.
12. Incubate the membrane in the diluted secondary antibody conjugate for 30 minutes at room temperature with shaking.
13. Wash the membrane with Wash Buffer four times for 5 minutes each.
14. Detect protein with the appropriate detection system.

**Table 1. Primary and secondary antibody concentrations to use with Thermo Scientific Chemiluminescent Substrates.**

Substrate	Pierce ECL	SuperSignal West Pico	SuperSignal West Dura	SuperSignal West Femto
<b>Primary Antibody Concentration (µg/mL)</b>	0.2-10	0.2-1.0	0.02-1.0	0.01-0.2
<b>Secondary Antibody Concentration (ng/mL)</b>	67-1000	10-50	4-20	2-10

## Troubleshooting

Problem	Possible Cause	Solution
Weak or no signal	Primary or secondary antibody concentration is too low	Use more primary or secondary antibody
	Too much enzyme in the system depleted the chemiluminescent substrate and caused the signal to fade quickly	Use less secondary antibody
	Inefficient protein transfer	Optimize transfer conditions
High background	Inadequate blocking and/or washing	Optimize blocking buffer and washing steps
	Secondary antibody conjugate concentration is too high	Use less secondary antibody conjugate

---

## Additional Information

Visit the website for additional information relating to this product including the following:

- Request your free copy of the Western Blotting Handbook, which contains a large troubleshooting section, detailed Western blotting protocols and dozens of SuperSignal Substrate references
- Tech Tip # 24: Optimize antigen and antibody concentrations for Western blots
- Tech Tip # 67: Chemiluminescent Western blotting technical guide and protocols
- Tech Tip # 21: Convert to SuperSignal West Pico Substrate from ECL

## Related Thermo Scientific Products

26681	Pierce Blue Prestained Protein Molecular Weight Marker Mix, 1 × 48 microtube plate
34080	SuperSignal West Pico Chemiluminescent Substrate, 500mL
34076	SuperSignal West Dura Extended Duration Substrate, 200mL
34096	SuperSignal West Femto Maximum Sensitivity Substrate, 200mL
32106	Pierce ECL Western Blotting Substrate, 500mL
37515	SuperBlock Blocking Buffer in PBS, 1L
37535	SuperBlock Blocking Buffer in TBS, 1L
21059	Restore™ Western Blot Stripping Buffer, 500mL
21065	Pierce Background Eliminator Kit, for eliminating background from X-ray film
34090	CL-XPosure™ Film, 5” × 7” sheets, 100 sheets/pkg

## General Reference

Alegria-Schaffer, A., *et al.* (2009). Performing and optimizing Western blots with an emphasis on chemiluminescent detection. *Methods Enzymol* **463**:573-99.

SuperSignal® Technology is protected by U.S. Patent # 6,432,662.

Tween® is a registered trademark of ICI Americas.

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.

© 2010 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.