

# SuperBlock<sup>®</sup> Dry Blend Blocking Buffer in TBS

37545

0705.4

Number	Description
37545	<b>SuperBlock Dry Blend Blocking Buffer in TBS</b> , 5 pouches, contains a proprietary protein formulation in Tris buffered saline, pH 7.4; each pouch yields 200mL

**Storage:** Upon receipt store at room temperature. After reconstitution, store at 4°C.

## Introduction

Thermo Scientific SuperBlock Blocking Buffer contains a proprietary non-relevant protein for blocking excess binding sites in ELISA, immunohistochemistry and Western blotting applications. This blocking buffer yields a high signal-to-noise ratio in most systems, does not contain albumin or endogenous biotin and, therefore, may be used with avidin/streptavidin systems.

## Important Product Information

- SuperBlock Blocking Buffer does not contain preservatives or anti-microbial agents. To prevent microbial contamination, anti-microbial agents may be added to the buffer or the buffer may be sterile filtered.
- Because no blocking reagent is optimal for all systems, empirical testing is essential to determine the appropriate blocking buffer for each Western blot system. Determining the proper blocking buffer can increase sensitivity and prevent nonspecific signal caused by cross-reactivity between the antibody and the blocking reagent.
- Reconstitute SuperBlock Blocking Buffer with 200mL of ultrapure water and use at the supplied concentration. However, other buffer concentrations may be beneficial for specific systems.
- A final concentration of 0.05% Tween<sup>®</sup>-20 Detergent added to the SuperBlock Blocking Buffer may improve performance; however, it is not required nor recommended for all systems. Use only high-quality products such as Thermo Scientific Surfact-Amps 20 (Product No. 28320), which is a specially purified Tween<sup>®</sup>-20 Detergent that is free of peroxides and carbonyls that may interfere in some systems.
- SuperBlock Blocking Buffer diluted 10-fold and containing 0.05% Tween<sup>®</sup>-20 Detergent may be used as a diluent for antibodies to improve signal-to-noise ratios.
- SuperBlock Blocking Buffer may be used as a protein stabilizer for drying antigen or antibody coated microplates. Dry plate completely before sealing in a plastic bag with desiccant. Store plate at 4°C.

## Procedure for Blocking ELISA Plates

1. Coat the ELISA plate with antigen or antibody.
2. Add 200µL of SuperBlock Blocking Buffer to each well.
3. Immediately empty the plate by inversion.
4. Repeat Steps 2-3 two additional times.
5. Proceed with assay or invert plate and allow it to completely dry. Place plate in a plastic bag or other container with desiccant and store at 4°C.

---

## Procedure for Blocking Membranes

1. Remove the membrane from the transfer unit.
2. Add enough SuperBlock Blocking Buffer to completely cover the membrane and incubate for 10-120 minutes with agitation.
3. Continue the Western blotting detection procedure.

## Procedure for Immunohistochemical Blocking

1. Block nonspecific sites in the tissue with SuperBlock Blocking Buffer for 30 minutes at room temperature.
2. Pour off the SuperBlock Blocking Buffer (do not rinse the tissue)
3. Continue the immunohistochemical detection procedure.

## Additional Information

Please visit the web site for additional information including the following items:

- Tech Tip #24: Optimize antigen and antibody concentrations for Western blots
- Tech Tip #67: Chemiluminescent Western blotting technical guide and protocols
- Tech Tip #32: Guide to enzyme substrates for Western blotting
- Tech Tip #33: Guide to enzyme substrates for ELISA

## Related Thermo Scientific Products

For a complete selection of quality reagents for immunochemical applications, please see our catalog or web site.

<b>25200-44</b>	<b>Precise™ Protein Gels</b> (see catalog or web site for a complete listing)
<b>34080</b>	<b>SuperSignal® West Pico Chemiluminescent Substrate</b> , 500mL
<b>34075</b>	<b>SuperSignal West Dura Extended Duration Substrate</b> , 100mL
<b>34095</b>	<b>SuperSignal West Femto Maximum Sensitivity Substrate</b> , 100mL
<b>21059</b>	<b>Restore™ Western Blot Stripping Buffer</b> , 500mL
<b>21065</b>	<b>Pierce® Background Eliminator Kit</b>
<b>34090</b>	<b>CL-XPosure™ Film</b> , 5" × 7" sheets, 100 sheets/pkg
<b>28376</b>	<b>BupH™ Tris Buffered Saline Packs</b> , 40 packs
<b>28372</b>	<b>BupH Phosphate Buffered Saline Packs</b> , 40 packs
<b>34028</b>	<b>1-Step Ultra TMB-ELISA</b> , 250mL
<b>37070</b>	<b>SuperSignal ELISA Pico Chemiluminescent Substrate</b> , 100mL
<b>34095</b>	<b>SuperSignal ELISA Femto Maximum Sensitivity Substrate</b> , 100mL

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.

© 2012 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.