

# SEA BLOCK Blocking Buffer

37527

0672.2

Number	Description
37527	<b>SEA BLOCK Blocking Buffer</b> , 500mL, contains steelhead salmon serum in phosphate-buffered saline and 0.1% sodium azide

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

## Introduction

The Thermo Scientific SEA BLOCK Blocking Buffer contains steelhead salmon serum, which does not interact significantly with mammalian antibodies and, therefore, results in minimal or no background. This buffer is ideal for blocking excess binding sites in ELISA, blotting, immunohistochemistry and other immunochemical applications.

## General Procedure

1. Add blocking solution to plate wells, blotting membrane or immunohistochemical slide.  
**Note:** If desired, the blocking buffer may be diluted with phosphate-buffered saline. For best results, empirically determine the optimal concentration to use.
2. Incubate 30 minutes to 2 hours at room temperature or 37°C.

## Related Thermo Scientific Products

We offer a wide range of products for immunochemistry. For a complete selection of high-quality reagents for most aspects of ELISA, Western blotting, immunohistochemistry and other immunotechnology applications see our catalog or web site.

Number	Description
15041	<b>Pierce® 96-Well EIA Plates - Corner Notch</b> , for ELISAs, 100 plates
15031	<b>Pierce 8-Well EIA Strip Plates - Clear, Corner Notch</b> , for ELISAs, 100 plates
15082	<b>MicroTube Racked System</b> , 10 racks of 96 individual 1.1mL tubes per rack
15075	<b>Reagent Reservoirs</b> , 50mL capacity, sterile, 200/pkg
28382	<b>BupH™ Carbonate-Bicarbonate Buffer Packs</b> , 40 packs
28372	<b>BupH Phosphate Buffered Saline Packs</b> , 40 packs
15100	<b>Pierce Maleic Anhydride Activated Polystyrene Strip Plates</b> , 5 plates
15110	<b>Pierce Maleic Anhydride Activated Polystyrene 96-Well Plates</b> , 5 plates

---

## Additional Information

Visit our web site for additional information relating to this product including the following:

- Tech Tip #67: Chemiluminescent Western blotting technical guide and protocols
- Tech Tip #21: Convert to SuperSignal<sup>®</sup> West Pico Substrate from ECL Substrate
- Tech Tip #23: Strip and reprobe Western blots
- Tech Tip #24: Optimize antigen and antibody concentrations for Western blots
- Tech Tip #32: Guide to enzyme substrates for Western blotting
- Tech Tip #43: Protein stability and storage
- Tech Tip #16: Block endogenous biotin
- Tech Tip #33: Guide to enzyme substrates for ELISA
- Western Blotting Handbook and Troubleshooting Guide

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.