

# Fish Serum Blocking Buffer

Catalog Number 37527

Pub. No. MAN0011315 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](https://www.thermofisher.com/support).

## Product description

Thermo Scientific™ Fish Serum Blocking Buffer (Cat. No. [37527](#)) is an ideal buffer to block excess binding sites in ELISA, western blotting, immunohistochemistry, and other immunochemical detection methods that involve mammalian samples. The buffer contains steelhead salmon serum, which lacks specific binding interactions with mammalian proteins, resulting in little to no background.

Fish Serum Blocking Buffer is supplied in a ready-to-use format. Depending on the system, the buffer can also be diluted with phosphate-buffered saline (PBS) if needed. For best results, empirically determine the optimal concentration before use.

## Contents and storage

Table 1 Fish Serum Blocking Buffer (Cat. No. [37527](#))

Contents	Amount	Storage
Fish Serum Blocking Buffer <sup>[1]</sup>	500 mL	2–8°C

<sup>[1]</sup> The buffer contains steelhead salmon serum in PBS and 0.1% sodium azide

## Block ELISA plates

1. Coat an ELISA plate with antigen or antibody.
2. Add 300 µL of Fish Serum Blocking Buffer to each well, then incubate for 30 minutes to 2 hours at room temperature or 37°C.
3. Remove the blocking buffer from the wells by aspiration or inversion.
4. Continue with the ELISA protocol that is appropriate for your downstream detection.

## Block immunohistochemistry tissue

1. Add an appropriate volume of Fish Serum Blocking Buffer to the tissue, then incubate for 30 minutes to 2 hours at room temperature or 37°C.
2. Pour off the blocking buffer. Do not rinse the tissue.
3. Continue with the immunohistochemical detection procedure that is appropriate for your downstream detection.

## Block western blots

1. After protein transfer, remove the membrane from the transfer apparatus, then wash in deionized water for 5 minutes, using agitation to remove all transfer buffer.
2. Add a sufficient amount of Fish Serum Blocking Buffer to cover the membrane.
3. Incubate for 30 minutes to 2 hours at room temperature with shaking.
4. Continue with the western blotting procedure that is appropriate for your downstream detection.

## Related products

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com).

Product	Cat. No.
ELISA reagents and kits	<a href="https://www.thermofisher.com/ELISA">thermofisher.com/ELISA</a>
ELISA plate readers	<a href="https://www.thermofisher.com/microplatereaders">thermofisher.com/microplatereaders</a>
Immunohistochemistry (IHC) reagents	<a href="https://www.thermofisher.com/cellimaging">thermofisher.com/cellimaging</a>
Western blotting reagents and accessories	<a href="https://www.thermofisher.com/westernblot">thermofisher.com/westernblot</a>

## 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](https://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](https://www.thermofisher.com/support).



**Revision history:** Pub. No. MAN0011315

Revision	Date	Description
B.0	20 July 2022	<ul style="list-style-type: none"><li>The user guide was updated to the current document template, with associated updates to the warranty, trademarks, and logos.</li><li>The product name was changed from SEA BLOCK Blocking Buffer to Fish Serum Blocking Buffer.</li></ul>
A.0	13 January 2012	Initial release.

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.

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

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.