

SuperBlock™ Blocking Buffers

Pub. No. MAN0011288

Rev. B.0

Pub Part No. 2160557.9

Number	Description
37580	SuperBlock Blocking Buffer in PBS , 100mL, contains a proprietary protein in phosphate-buffered saline (pH 7.4) containing Kathon™ Anti-microbial Agent
37515	SuperBlock Blocking Buffer in PBS , 1L, contains a proprietary protein in phosphate-buffered saline (pH 7.4) containing Kathon Anti-microbial Agent
37516	SuperBlock T20 (PBS) Blocking Buffer , 1L, contains a proprietary protein formulation in phosphate buffered saline (pH 7.4) containing 0.05% Tween™-20 Detergent and Kathon Anti-microbial Agent
37518	SuperBlock Blocking Buffer in PBS , 5L, contains a proprietary protein in phosphate-buffered saline (pH 7.4) containing Kathon Anti-microbial Agent
37581	SuperBlock Blocking Buffer in TBS , 100mL, contains a proprietary protein in Tris-buffered saline (pH 7.4) containing Kathon Anti-microbial Agent
37535	SuperBlock Blocking Buffer in TBS , 1L, contains a proprietary protein in Tris-buffered saline (pH 7.4) containing Kathon Anti-microbial Agent
37536	SuperBlock T20 (TBS) Blocking Buffer , 1L, contains a proprietary protein formulation in Tris-buffered saline (pH 7.4) containing 0.05% Tween-20 Detergent and Kathon Anti-microbial Agent

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

Introduction

The Thermo Scientific SuperBlock Blocking Buffers contain 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 and does not contain albumin or endogenous biotin and, therefore, may be used with avidin/streptavidin systems. For ease of use, SuperBlock T20 Blocking Buffers contain Tween-20 Detergent, which improves blocking performance in many Western blot detection systems.

Important Product Information

- 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.
- Use SuperBlock Blocking Buffers as supplied. However, other buffer concentrations may be beneficial for specific systems. For example, when using the blocking buffers as a diluent for antibodies to improve signal-to-noise ratios, the buffers may be used as supplied or diluted up to 10-fold.
- A final concentration of 0.05% Tween-20 in the blocking buffer often improves blocking; however, it is not required nor recommended for all systems. Use only high-quality products such as Thermo Scientific Surfact-Amps™ 20 Detergent Solution (Product No. 28320), which is a specially purified Tween-20 Detergent that is free of peroxides and carbonyls that may interfere in some systems. SuperBlock T20 Blocking Buffers are supplied containing Tween-20 Detergent at a concentration of 0.05%.
- SuperBlock Blocking Buffers 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 300µL of blocking buffer to each well and immediately empty the plate. Repeat this step two additional times.
3. Proceed with assay or invert plate and allow it to completely dry for ~2 hours. 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 blocking buffer to cover the membrane and incubate for 10-120 minutes with agitation.
3. Continue the Western blot detection procedure. Use the blocking buffer to dilute primary and secondary antibodies.

Procedure for Blocking Tissue for Immunohistochemistry

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

Additional Information from Our Website

- Tech Tip #24: Optimize antigen and antibody concentrations for Western blots
- Tech Tip #16: Block endogenous biotin

Related Products

XP04200BOX	Novex™ Tris-Glycine protein gels (see thermofisher.com/proteingels for a complete listing)
NW04120BOX	Bolt™ Bis-Tris Plus protein gels (see thermofisher.com/proteingels for a complete listing)
28376	BupH™ Tris Buffered Saline Packs, 40 packs, each pack yields 500mL
28372	BupH Phosphate Buffered Saline Packs, 40 packs, each pack yields 500mL
34580	SuperSignal™ West Pico Plus Chemiluminescent Substrate, 500mL
34075	SuperSignal West Dura Extended Duration Substrate, 100mL
34095	SuperSignal West Femto Maximum Sensitivity Substrate, 100mL
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
37527	SEA BLOCK Blocking Buffer, 500mL
37582	Blocker Casein in PBS, 100 mL
37528	Blocker Casein in PBS, 1L
37583	Blocker Casein in TBS, 100 mL
37530	Blocker BLOTTO in TBS, 1L
37584	Pierce Protein-Free (PBS) Blocking Buffer, 100 mL
37572	Pierce Protein-Free (PBS) Blocking Buffer, 1L
37573	Pierce Protein-Free T20 (PBS) Blocking Buffer, 1L
37585	Pierce Protein-Free (TBS) Blocking Buffer, 100 mL
37570	Pierce Protein-Free (TBS) Blocking Buffer, 1 L
37571	Pierce Protein-Free T20 (TBS) Blocking Buffer, 1 L

Tween is a trademark of Croda International Plc.

Kathon is a trademark of The Dow Chemical Company.

For research use only. Not for use in diagnostic procedures.

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 thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 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.