

StartingBlock™ Blocking Buffers

1457.4

Number	Description
37578	StartingBlock (PBS) Blocking Buffer , 100 mL, contains a proprietary protein formulation in phosphate-buffered saline at pH 7.5
37538	StartingBlock (PBS) Blocking Buffer , 1L, contains a proprietary protein formulation in phosphate-buffered saline at pH 7.5
37539	StartingBlock T20 (PBS) Blocking Buffer , 1L, contains a proprietary protein formulation in phosphate-buffered saline at pH 7.5 with 0.05% Tween®-20 Detergent
37579	StartingBlock (TBS) Blocking Buffer , 100 mL, contains a proprietary protein formulation in Tris-buffered saline at pH 7.5
37542	StartingBlock (TBS) Blocking Buffer , 1L, contains a proprietary protein formulation in Tris-buffered saline at pH 7.5
37543	StartingBlock T20 (TBS) Blocking Buffer , 1L, contains a proprietary protein formulation in Tris-buffered saline at pH 7.5 with 0.05% Tween-20 Detergent

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

Introduction

The Thermo Scientific StartingBlock Blocking Buffer is the first blocking buffer to use for optimizing a new system or when background signal is occurring in the current system. Robustness and lack of cross-reactivity are key characteristics for determining the optimal protein-based blocking buffer to eliminate binding sites in ELISA, blotting, immunohistochemistry and other immunochemical applications. StartingBlock Blocking Buffers are compatible with antibodies and biotin/avidin systems and results in high signal-to-noise ratios. For ease of use, StartingBlock T20 Blocking Buffer contains the detergent Tween-20, which improves blocking performance in many Western blot detection systems.

Important Product Information

- The usage as described in these instructions may differ from other blocking solutions.
- Use StartingBlock Buffers at the supplied concentration; do not dilute blocking buffer.

Procedure for Blocking Membranes for Western Blotting

Note: For best results, use StartingBlock T20 Blocking Buffer or add a final concentration of 0.05% Tween-20 Detergent to StartingBlock Blocking Buffer.

1. Remove the membrane from the transfer unit. Add sufficient blocking buffer to cover the membrane.
2. For direct methods (i.e., using an enzyme-labeled primary antibody), incubate blot for 1-5 minutes at room temperature with agitation. For indirect methods (i.e., using an enzyme-labeled secondary), incubate for 15-30 minutes at room temperature with agitation.
3. Continue the Western blotting procedure.

Note: For best results, use StartingBlock T20 Blocking Buffer or StartingBlock Blocking Buffer containing a final concentration of 0.05% Tween-20 Detergent as an antibody diluent. If the blocker is not used as a diluent for direct methods, blocking time may need to be increased to 10-15 minutes.

Procedure for Blocking Microplates for ELISA

Note: Generally, it is not necessary to add Tween-20 to the blocking buffer when used for blocking. However, using StartingBlock T20 Blocking Buffer when used as an antibody diluent may increase the signal-to-noise ratio.

1. Coat the ELISA plate with antigen or antibody.
2. Add 300 μ L of blocking buffer to each well. Empty plate by aspiration or inversion. Incubation is not required before emptying plate.
3. Proceed with the ELISA protocol or, for storage, invert plate for approximately 2 hours to dry.
4. Transfer plate to a plastic bag or other container containing a desiccant, such as silica gel. Store the plate at 4°C.

Note: Most antibodies and antigens are stable for greater than one year when coated onto plates and stored dry.

Additional Information from Our Website

- Tech Tip #24: Optimize antigen and antibody concentrations for Western blots
- Tech Tip #67: Chemiluminescent Western blotting technical guide and protocols
- Protein Methods Library: Blocking Buffers for Western Blotting and ELISA

Related Thermo Scientific Products

34080	SuperSignal[®] West Pico Chemiluminescent Substrate, 500mL
34075	SuperSignal West Dura Extended Duration Substrate, 100mL
34095	SuperSignal West Femto Maximum Sensitivity Substrate, 100mL
46430	Restore Plus Western Blot Stripping Buffer, 500mL
21059	Restore[™] Western Blot Stripping Buffer, 500mL
34090	CL-Xposure[™] Film, 5" × 7" sheets, 100 sheets/pkg
21065	Pierce[®] Background Eliminator Kit, for eliminating background from X-ray film
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
37580	SuperBlock[®] Blocking Buffer in PBS, 100 mL
37515	SuperBlock Blocking Buffer in PBS, 1L
37516	SuperBlock T20 (PBS) Blocking Buffer, 1L
37581	SuperBlock Blocking Buffer in TBS, 100 mL
37535	SuperBlock Blocking Buffer in TBS, 1L
37536	SuperBlock T20 (TBS) Blocking Buffer, 1L
37584	Pierce Protein-Free (PBS) Blocking Buffer, 100 mL
37585	Pierce Protein-Free (TBS) Blocking Buffer, 100 mL
28376	BupH[™] Tris Buffered Saline Packs, 40 packs
28372	BupH Phosphate Buffered Saline Packs, 40 packs
25200-74	Precise[™] Protein Gels

General Reference

Antibodies: A Laboratory Manual. Ed Harlow and David Lane, Cold Springs Harbor Laboratory, 1998.

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. 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.