

Blocker BSA (10X)

37520 37525

0551.4

Number	Description
37520	Blocker BSA in TBS (10X) , 125mL, contains 10% bovine serum albumin in 10mM Tris, 150mM NaCl, containing Kathon™ Anti-microbial Agent; pH 7.4
37525	Blocker BSA in PBS (10X) , 200mL, contains 10% bovine serum albumin in 10mM sodium phosphate, 150mM NaCl, containing Kathon Anti-microbial Agent; pH 7.4

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

Introduction

Proteinaceous solutions are used for blocking excess binding sites in ELISA, blotting, immunohistochemistry and other immunochemical applications. Each blocking solution will impart certain characteristics on the assay system. Robustness and lack of cross-reactivity are key characteristics in buffer selection. Thermo Scientific™ Blocker™ BSA is compatible with biotin/avidin systems and produces high signal-to-noise ratios in most systems due to its lack of endogenous biotin.

Important Product Information

- Because no blocking reagent is optimal for all systems, empirical testing is essential to determine the appropriate blocking buffer for each system. Determining the proper blocking buffer can increase sensitivity and prevent nonspecific signal caused by cross-reactivity between the antibody and the blocking reagent.
- The Blocker BSA (10X) is supplied as a 10% solution and should be diluted before use. Dilute using either 1X TBS for Product No. 37520 or 1X PBS for Product No. 37525. Solutions containing 1-3% BSA are used for most applications. Other BSA concentrations may be beneficial for specific systems.
- A final concentration of 0.05% Tween™-20 Detergent added to the Blocker BSA may improve blocking; 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-20 Detergent free of peroxides and carbonyls that may interfere in some systems.
- The Blocker BSA diluted 10-fold and containing 0.05% Tween-20 Detergent may be used as a diluent for antibodies to improve signal-to-noise ratios.
- Blocker BSA may be used as a protein stabilizer for drying antigen or antibody coated microplates. Perform the example procedure below and dry plate completely before sealing in a plastic bag with desiccant. Store plate at 4°C.

Example Procedure for Blocking

1. Dilute Blocker BSA to the appropriate concentration using 1X TBS or 1X PBS (see Important Product Information).
2. Add the blocking solution to the microplate well, blotting membrane or immunohistochemical slide.
3. Incubate 30 minutes to 2 hours at room temperature or 37°C.
4. Continue protocol for the specific application.

Additional Information on Our Website

- Tech Tip: Optimize antigen and antibody concentrations for Western blots
- Tech Tip: Determine source of nonspecific signal in chemiluminescent Western blots
- Tech Tip: Substrates for blotting and ELISA applications
- Tech Tip: Block endogenous biotin
- Tech Tip: Convert to SuperSignal™ West Pico Substrate from ECL Substrate

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
28376	BupH™ Tris Buffered Saline Packs, 40 packs
28372	BupH™ Phosphate Buffered Saline Packs, 40 packs
25200-44	Precise™ Protein Gels
88018	Nitrocellulose Membrane, 0.45µm, 30cm × 3.5m, roll
88024	Nitrocellulose Membrane, 0.2µm, 8cm × 8cm, 15/pkg
88025	Nitrocellulose Membrane, 0.45µm, 8cm × 8.5cm, 15/pkg
77012	Nitrocellulose Membrane, 0.2µm, 8cm × 12cm, 25/pkg
77010	Nitrocellulose Membrane, 0.45µm, 8cm × 12cm, 25/pkg
88585	PVDF (polyvinylidene difluoride) Membrane, 0.45µm, 10cm × 10cm, 10/pkg
88518	PVDF (polyvinylidene difluoride) Membrane, 0.45µm, 26.5cm × 3.75m, roll

Products are 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”). 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 Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

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 REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER’S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR 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.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com/techresources.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2015 Thermo Fisher Scientific Inc. All rights reserved. Kathon is a trademark of The Dow Chemical Company. Tween is a trademark of Croda International Plc. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.