Blocker[™] FL Fluorescent Blocking Buffer (10X)

Catalog Number 37565

Doc. Part No. 2160627 Pub. No. MAN0017084 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**.

Product description

Thermo Scientific[™] Blocker[™] FL Fluorescent Blocking Buffer (10X) is a single purified protein for blocking excess nonspecific binding sites to reduce background fluorescence in western blotting applications when used with nitrocellulose and PVDF membranes. This blocking buffer is ideal for fluorescent protein detection by providing lower background in blotting applications where background fluorescence may otherwise interfere with signal detection. The blocking solution is also compatible with chemiluminescent detection. Blocker[™] FL Fluorescent Blocking Buffer (10X) can be used for blocking of cells or tissue in fluorescence imaging.

Contents and storage

Product	Cat. No.	Amount	Storage
Blocker™ FL Fluorescent Blocking Buffer (10X)	37565	100 mL	4°C

Procedural guidelines

 Blocker[™] FL Blocking Buffer may be used at higher or more dilute concentrations. For best results, empirically determine the optimal concentration to use.

Block western blots

- 1. Prepare blocking buffer working reagent by adding 1 part Blocker[™] FL Fluorescent Blocking Buffer (10X) to 9 parts ultrapure water. Prepare sufficient volume of working reagent to completely cover the membrane.
- 2. After the protein transfer, remove the membrane from the transfer apparatus, then wash in deionized water for 5 minutes, using agitation to remove all transfer buffer.
- 3. Add sufficient Blocker[™] FL Fluorescent working reagent to cover the membrane.
- 4. Incubate 15-30 minutes at room temperature with shaking.
- 5. Continue with the western blotting procedure that is appropriate for your downstream detection.



Related products

Table 1 Blocking buffers for western blotting

Select when	Product	Blocking agent	Highlights	Available formats
Optimizing a new western blot system or high background with current blocking buffer	StartingBlock™ Blocking Buffer	Serum- and biotin-free, single purified protein	Performs well with a wide range of antibodies and antibody combinations Compatible with streptavidin systems Blocks in less than 15 minutes	PBS TBS PBST TBST
Use when high background or antigen-antibody masking occurring with non-fat milk blockers	Blocker [™] Casein	Purified casein	Single purified protein blocking buffer provides fewer chances of cross- reaction with assay components than serum or milk solutions	PBS TBS
Fluorescent western blotting	Blocker™ FL Fluorescent Blocking Buffer	Single purified protein	Blocks excess nonspecific sites to help reduce background fluorescence Detergent-free Blocks in 15-30 minutes	10X concentrate

Table 2 Additional products

Products	Learn more
Western blotting reagents and accessories	thermofisher.com/westernblot
Western blot imaging and analysis	thermofisher.com/westernimaging

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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.

Revision history: Pub. No. MAN0017084

Revision	Date	Description
B.0	17 September 2021	Updated format
A.0	20 June 2017	New document

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

©2021 Thermo Fisher Scientific Inc. All rights reserved. Kathon is a trademark of DuPont Chemical Company. Tween is a trademark of Croda International Plc. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

