

# Blocker FL 10X Fluorescent Blocking Buffer

MAN0017084

Rev. A.0

Pub. Part No. 2160672.3

## 37565

Number	Description
37565	Blocker FL 10X Fluorescent Blocking Buffer, 100mL

**Storage:** Upon receipt store at 4°C. Product is shipped on ice pack.

**For Research Use Only. Not for use in diagnostic procedures.**

## Introduction

The Thermo Scientific™ Blocker™ FL 10X Fluorescent Blocking Buffer is a proprietary formulation 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 10X Fluorescent Blocking Buffer can be used for blocking of cells or tissue in fluorescence imaging.

## General Procedure

1. Prepare working reagent by adding 1 part Blocker FL 10X to 9 parts ultrapure water. Prepare sufficient volume of working reagent to completely cover the membrane (recommended volume of 1mL working reagent per 10cm<sup>2</sup> of membrane).

**Note:** If desired, the blocking buffer may be used at a higher or more dilute concentrations. For best results, empirically determine the optimal concentration to use.

2. Add blocking solution working reagent to blotting membrane.
3. Incubate 15 to 30 minutes at room temperature.

## Related Thermo Scientific Products

We offer a wide range of products for Western blotting. For a complete selection of high-quality reagents for most aspects of ELISA, Western blotting, immunohistochemistry and other immunotechnology applications see our catalog or website.

Catalog no.	Name	Size
A32752	iBright™ FL1000 Instrument	1 unit
A32749	iBright™ CL1000 Instrument	1 unit
LC5615	iBright™ Prestained Protein Ladder	2 × 250µL
LC2570	Fluorescent Compatible Sample Buffer (4X), Non-Reducing	5mL
NW0412A	Bolt™ Welcome Pack A (4-12%, 10-well)	1 kit
NW0412B	Bolt Welcome Pack A (4-12%, 15-well)	1 kit
22860	Low Fluorescence PVDF Transfer Membrane, 7 cm x 8.4 cm	10 sheets
A32723	Goat anti-Mouse IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor™ Plus 488	1mg
A32727	Goat anti-Mouse IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 555	1mg
A32728	Goat anti-Mouse IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 647	1mg
A32729	Goat anti-Mouse IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 680	1mg
A32730	Goat anti-Mouse IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 800	1mg
A32731	Goat anti-Rabbit IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 488	1mg
A32732	Goat anti-Rabbit IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 555	1mg
A32733	Goat anti-Rabbit IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 647	1mg
A32734	Goat anti-Rabbit IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 680	1mg
A32735	Goat anti-Rabbit IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa Fluor Plus 800	1mg

---

## Additional Information

Visit our website for additional information relating to this product including the following:

- Tech Tip #67: Chemiluminescent Western blotting technical guide and protocols
- Tech Tip #24: Optimize antigen and antibody concentrations for Western blots
- Tech Tip #32: Guide to enzyme substrates for Western blotting
- Tech Tip #43: Protein stability and storage
- Tech Tip #16: Block endogenous biotin
- Western Blotting Handbook and Troubleshooting Guide

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](http://thermofisher.com). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.