



043PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# BLOT-QuickBlocker™

(Cat. # 786-011)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Protein solutions are used for blocking excess binding sites in western blots and other immunochemical applications. Each blocking solution will impart certain characteristics on the assay system. BLOT-QuickBlocker™ is a modified milk protein, and is good for blocking western blot membranes.

## ITEM(S) SUPPLIED (Cat. # 786-011)

| Description        | Size |
|--------------------|------|
| BLOT-QuickBlocker™ | 175g |

## STORAGE CONDITIONS

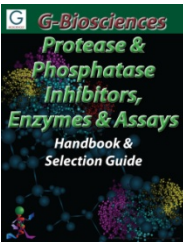
It is shipped at ambient temperature. Store at room temperature, stable for 1 year.

## PROTOCOL

1. Prepare 5% solution of BLOT-QuickBlocker™ in TBS /TBST or PBS/PBST.
2. Add 15-20ml of 5% solution of BLOT-QuickBlocker™ on to the blotting membrane. Make sure the membrane is submerged in the solution.
3. Incubate the membrane with BLOT-QuickBlocker™ for 30-60 minutes or more at room temperature with gentle shaking.
4. Proceed with the blot detection protocol as per your specific application.

## RELATED PRODUCTS

Download our Protease & Phosphatase Inhibitors, Enzyme & Assays Handbook.



<http://info.gbiosciences.com/protease-phosphatase-inhibitors-enzymes-assay-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Last saved: 7/30/2012 CMH

*This page is intentionally left blank*



[www.GBiosciences.com](http://www.GBiosciences.com)