



035PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

BLOK™-BLOTTO in TBS

(Cat. # 786-192)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

BLOK™-BLOTTO is a pre-formulated, non-fat milk solution for blocking excess binding sites in ELISA, Western blotting, immunohistochemistry and other immunochemical applications. It contains 5% (w/v) nonfat powdered milk in 25mM Tris, 150mM NaCl, pH 7.4 with 0.01% Antifoam A in ready to use form.

ITEM(S) SUPPLIED (Cat. # 786-192)

Description	Size
BLOK™-BLOTTO in TBS (1X)	2 x 500ml

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store it refrigerated at 4°C and is stable for 1 year.

PROTOCOL

1. Gently swirl the bottle.
2. Add the BLOK™-BLOTTO in TBS solution to the microplate well, blotting membrane or immunohistochemical slide.
3. Incubate 0.5-2 hours at room temperature with gently shaking.
4. Continue with protocol for your specific application.

RELATED PRODUCTS

Download our Western Blotting and Assay Development Handbooks



<http://info.gbiosciences.com/complete-western-blot-handbook--selection-guide>

<http://info.gbiosciences.com/complete-assay-development-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 11/12/2012 CMH

This page is intentionally left blank



www.GBiosciences.com