



039PR

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

A Geno Technology, Inc. (USA) brand name

BLOK™ -BSA in TBS

(Cat. # 786-193)



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

INTRODUCTION

Protein 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. BLOK™-BSA in TBS does not contain endogenous biotin, therefore it is compatible with biotin/avidin systems and results in high signal-to-noise ratios in most systems. It is robust and shows a lack of cross-reactivity.

ITEM(S) SUPPLIED (Cat # 786-193)

Description	Size
BLOK™-BSA in TBS [10X]	125ml

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store at 4°C. When stored and used properly, the kit is stable for 1 year.

PROTOCOL

1. Gently swirl the bottle. Dilute appropriate volume of the BLOK™-BSA in TBS 10 fold with 1X TBS (e.g. mix 2ml BLOK™-BSA with 18ml 1X TBS). The final concentration of BSA will be 1%.
2. Add the 1X BLOK™-BSA in TBS solution into the micro-plate well, blotting membrane or immunohistochemical slide.
3. Incubate for 30 minutes to 2 hours at room temperature with gentle shaking.
4. Proceed to the 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.

This page is intentionally left blank



www.GBiosciences.com