



037PR-02

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

A Geno Technology, Inc. (USA) brand name

BLOK™ BSA in PBS

(Cat. # 786-195)



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 PBS 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. # 786-195)

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

STORAGE CONDITIONS

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

PROTOCOL

1. Gently swirling the bottle. Dilute appropriate volume of the BLOK™ BSA in PBS 10 fold with 1X PBS (e.g. mix 2ml BLOK™ BSA in PBS with 18ml 1X PBS). The final concentration of BSA will be 1%.
2. Add the 1X BLOK™ BSA in PBS solution to 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