



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

A Geno Technology, Inc. (USA) brand name

Superior™ Blocking Buffer for Precipitating Substrate

For Blocking Application with Precipitating
Substrates on Membranes

(Cat. # 786-655, 786-656)



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

INTRODUCTION

Superior™ Blocking Buffer for Precipitating Substrate contains a proprietary antigenically non-determinant protein for blocking non-specific sites during ELISA, membrane blotting, immunohistochemistry and other applications. This blotting buffer has been optimized for used in blotting protocols that use precipitating substrates. Superior™ Blocking Buffer is ideal for a high signal to background ratio in most system. Superior™ Blocking Buffer does not contain biotin or other animal source proteins to interfere with immuno-complexes. Superior™ Blocking Buffer is suitable for assays that use avidin/streptavidin systems.

ITEM (S) SUPPLIED

Cat. #	Description	Size
786-655	Superior™ Blocking Buffer for Precipitating Substrate in TBS (Tris-buffered saline at pH 7.5)	500ml
786-656	Superior™ Blocking Buffer for Precipitating Substrate in PBS (Phosphate-buffered saline at pH 7.5)	500ml

STORAGE CONDITIONS

Shipped at ambient conditions, upon arrival store at 4°C.

IMPORTANT

- For optimal blocking, do NOT dilute the Superior™ Blocking Buffer.
- This product is specifically formulated for precipitate enzyme dye substrate and it is not suitable for chemiluminescent or ELISA application. See related product for other applications.
- The efficacy of blocking agents varies from application to application, so we recommend empirical testing of blocking buffer and optimization of procedure to increase sensitivity and prevent nonspecific signal and cross-reaction between blocking agent and antibody.
- Use of detergent in blocking buffers is not required for all applications, however, addition of 0.05% Tween®-20 often improves blocking. Use only high quality ultra pure grade Tween®-20, we recommend our Proteomic Grade Tween®-20 solution (Cat. # DG011, DG012, DG511), which is purified to remove peroxide and carbonyl contaminants that may interfere in some applications.
- 10-fold diluted Superior™ Blocking Buffer containing 0.05% Tween®-20 may be used to dilute antibodies to enhance the sensitivity of the signal.

PROCEDURE FOR BLOCKING WESTERN BLOTTING MEMBRANES

1. Following protein transfer to the membrane, transfer the membrane to a suitable size tray.

NOTE: Superior™ Blocking Buffer is suitable for PVDF and nitrocellulose membranes.

2. Add enough Superior™ Blocking Buffer to completely cover the membrane.
3. Incubate for 30-120 minutes at room temperature with agitation.
4. Discard blocking buffer and continue with downstream Western blotting steps.

NOTE: For washing steps, use of femtoTBST™ (Cat. # 786-161) or femtoPBST™ (Cat. # 786-162) will minimize the washing out of immune-complexes and aid in the generation of cleaner backgrounds resulting in a higher signal to noise ratio, a common problem associated with classical TBST and PBST buffers used for washing.

RELATED PRODUCTS

Download our Western Blotting Handbook and Protein Discovery.



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

<http://info.gbiosciences.com/complete-protein-discovery-handbook>

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

Last saved: 8/15/2013 SM



www.GBiosciences.com