



041PR-01

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

A Geno Technology, Inc. (USA) brand name

BLOT-FastStain™

Rapid, Sensitive and Reversible Staining
of Proteins on Western Blots

(Cat. # 786-34)



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

INTRODUCTION

BLOT-FastStain™ is a unique proprietary system for the reversible staining of proteins after transfer to nitrocellulose or PVDF membranes. BLOT-FastStain™ detects only proteins, and leaves the background absolutely untouched and brilliant white, leading to an exceptional band visibility. The lower detection limit of BLOT-FastStain™ is 0.5ng protein (BSA)/band on nitrocellulose membrane.

Perhaps more importantly than simply detecting proteins on transfer membranes, BLOT-FastStain™ can be removed in 10 minutes without affecting their biological or immunological properties. After removing the stain, Western blotting can proceed with no changes in protocol. BLOT-FastStain™, therefore, allows you to measure the transfer of proteins before potentially wasting expensive and sometimes irreplaceable antibodies. Other applications include detecting size-fractionated proteins for later antibody generation or sequencing.

ITEM(S) SUPPLIED (Cat. # 786-34)

Description	Size
Reagent A: Fixer	1 x 120 ml
Reagent B: Developer	2 x 125 ml

The kit components are sufficient for 25 blots of 12cm x 12cm size.

STORAGE & STABILITY

The kit is shipped at ambient temperature. Upon arrival, store reagents at room temperature. Do not store in the cold. The kit components are stable for 12 months, when stored and handled properly. To avoid possible microbial contamination, dispense all solutions aseptically.

STAINING PROTOCOL

Protocol is for a single 12 x 12cm membrane, adjust volumes proportionally for larger membranes.

1. Dilute Reagent A: Fixer 10 fold by adding 4ml of the fixer to 36 ml DI water. Place transfer membrane (nitrocellulose or PVDF) into a tray containing 40ml diluted fixer. Incubate for 2-3 minutes with gentle shaking.
2. Dilute Reagent B: Developer 4 fold by adding 10ml of the supplied Developer to 30 ml DI water. Transfer the membrane (nitrocellulose or PVDF) into the tray containing 40ml diluted Developer. Incubate for 1 minute with gentle shaking.
3. Transfer the Developer tray into a refrigerator for protein bands to develop. Shaking at this stage is not recommended.
4. Incubate until protein bands can be detected and reach their desired intensity. The best results will be achieved if the membranes are developed in the cold and dark condition without shaking or disturbing.

Protein bands will be visible within 10 minutes of incubation and reaches full intensity within 30 minutes. Developing time depends on the concentration of protein on the surface of the membrane. After staining, membrane may be left in the developer for several days.

BACKGROUND STAINING

Background staining depends on the types of membrane used for protein transfer. Nitrocellulose gives the clearest and most brilliant white background. Some PVDF membranes may give higher background. Background staining can be removed by shaking the membrane 5-10 minutes in cold water (too much washing may de-stain the protein bands).

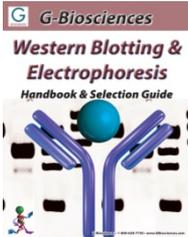
DESTAINING

To destain, simply place the membrane in warm deionized water (40-45°C) and shake in bright light for 10 minutes or until stain disappears. Wash twice, 10 minute each, with deionized water.

Most bands will fade away within 10 minutes. Some protein bands may take longer, depending on their protein concentration as well as their amino acid composition.

RELATED PRODUCTS

Download our Western Blotting Handbook.



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

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

Last saved: 10/3/2012 CMH



www.GBiosciences.com