



140PR

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

A Geno Technology, Inc. (USA) brand name

Coomassie Brilliant Blue R-250 Stain

For Staining of Polyacrylamide Gels

(Cat. # 786-498)



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

INTRODUCTION

Coomassie Brilliant Blue R-250 Stain is based on a colloidal Coomassie stain. Coomassie Brilliant Blue R-250 Stain stains proteins, with high band visibility. The staining of gels with Coomassie Brilliant Blue R-250 Stain allows the examination of protein bands even during the staining process. After the staining process, the band intensity may be further enhanced by de-staining the stained gel in our Coomassie Brilliant Blue Destaining Solution (Cat# 786-499) or 30% Methanol. Coomassie Brilliant Blue R-250 Stain has a sensitivity of staining 8-10ng protein/band, i.e. 8-10ng of BSA is visible in 4-20% SDS acrylamide gels. The supplied volume is suitable for 40-50 Mini gels (8 x 10cm) & 20-25 large gels (12 x 15cm).

ITEM(S) SUPPLIED (CAT. # 786-498)

Description	Size
Coomassie Brilliant Blue R-250 Stain	1L

STORAGE CONDITION

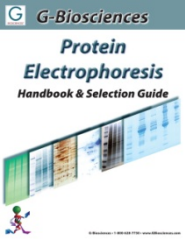
It is shipped at ambient temperature. Store it at room temperature, upon arrival. It is stable for 1 year when stored and used properly.

PROTOCOL FOR STAINING THE POLYACRYLAMIDE GELS:

1. Wash gel 2-3 times, 5 minutes each in deionized water to remove SDS present in the gel. Each wash should be in large volumes of water. For gels without SDS, a single wash in deionized water is sufficient.
NOTE: *Isoelectric focusing gels require prefixing in 20% trichloroacetic acid for 30 minutes, followed by extensive washing to remove TCA.*
2. Remove all free water from the gel. Add an adequate amount of Coomassie Brilliant Blue R-250 stain to cover the gel. Gently shake the gel in stain for 1 hour. Protein bands will be visible within 3-5 minutes and reach a maximum intensity within 1 hour. Longer incubation may be performed.
3. Rinse the stained gel in a large volume of deionized water, 2-3 times for 5 minutes each. Rinsing in deionized water enhances the intensity of the protein bands. Rinse gel in Coomassie Brilliant Blue De-Staining Solution (Catalog # 786-499) or 30% Methanol, until desired resolution is attained. Store the stained gel in deionized water.
NOTE: *If background staining is noticed, it is indicative of residual SDS in the gel. Rinsing the gel extensively in deionized water will remove the background staining.*

RELATED PRODUCTS

Download our Protein Electrophoresis Handbook



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

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

Last saved: 8/16/2012 CMH



www.GBiosciences.com