



381PR

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

A Geno Technology, Inc. (USA) brand name

Silver D-Stain™

(Cat. # 786-199)



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

INTRODUCTION

Silver staining is one of the most sensitive methods for detecting proteins and nucleic acids. One of the most common problems associated with silver staining of gels is detection and resolution of both weak as well as strong bands and spots on the gel. If a gel is developed for detecting stronger bands or spots, the weaker bands or spots are either hard to detect or not detected at all. If, on the other hand, a gel is developed to detect weaker bands or spots, the stronger bands or spots will become too dark, obscuring resolution of the entire gel. The Silver D-Stain™ is specifically developed for detecting all of the images in any gel - weaker as well as stronger signals. The Silver D-Stain™ will allow either complete or partial destaining of acrylamide gels. The destained or partially destained gels are ready for any downstream applications, including re-staining with any method to obtain satisfactory resolution of the gel. The Silver D-Stain™ is suitable for both protein and nucleic acid acrylamide gels. The Silver D-Stain™ kit is sufficient for processing at least 25 mini gels.

ITEM(S) SUPPLIED (Cat. # 786-199)

Description	Size
Silver D-Stain™ Part-I	250ml
Silver D-Stain™ Part-II	250ml

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store at room temperature. When stored and handled properly this kit is stable for 1 year.

PROTOCOL

1. Wash the silver stained gel 3-4 times, 10 minutes each, with generous amount of de-ionized water.
2. Prepare a working reagent solution - mix equal volumes of Part-I and Part-II.
NOTE: Prepare sufficient volume of reagent solution to completely submerge the gel to be destained. For a mini gel, mix 10 ml from each part. For larger gels, increase the volume as necessary.
For partially destaining of a gel or for allowing the destaining to proceed at a slower rate for better control, dilute the working reagent solution by 3-5 fold with pure water.
3. Transfer the gel into working reagent. Gently rock the gel.
4. Observe the gel for destaining. Destaining may be stopped at any stage or when desired result is achieved. Destaining is generally complete within 5 - 10 minutes. Some strong bands/spots may take longer to destain.
5. When the gel has reached the desired resolution, quickly rinse the gel twice with de-ionized water. Wash the gel three times, 5-10 minutes each with water. The gel is now ready for next step.

RESTAINING GELS.

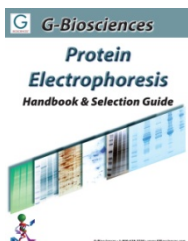
The destained gel is ready for downstream processing, including restaining, either with silver stain (*FastSilver™*, Cat. # 786-30) or any other method.

PREPARATION FOR WESTERN BLOTTING

If the destained gel is to be used for transfer of protein or nucleic acids to transfer membranes (Western or Northern Transfer), before transfer soak the gel in generous amount of transfer buffer for 30-50 minutes.

RELATED PRODUCTS

Download our Protein Cross-linkers 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: 2/28/2013 CMH



www.GBiosciences.com