



580PR-01

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

A Geno Technology, Inc. (USA) brand name

SilverOUT™

**For the Removal of Silver Ions
Following Silver Staining**

(Cat. # 786-244)



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

INTRODUCTION

SilverOUT™ is a two component system designed for the removal of silver ions for protein gel species stained with a silver protein stain (G-Biosciences, FASTsilver™ (Cat. # 786-30) or FOCUS™ FASTsilver™ (Cat. # 786-240).

The destained gel pieces are suitable for further downstream applications, including reduction, alkylation and peptide digestion with trypsin for mass spec analysis.

The kit is suitable for >100 protein gel spots.

ITEM(S) SUPPLIED (Cat. # 786-244)

Description	Size
SilverOUT™ I	4ml
SilverOUT™ II	4ml

STORAGE CONDITION

Shipped at ambient temperature. Store at 4°C upon arrival.

ADDITIONAL ITEMS REQUIRED

- Ultrapure water (18MΩ equivalent), we recommend our Proteomic Grade Water (Cat. # 786-229)

IMPORTANT INFORMATION

To reduce keratin and chemical backgrounds

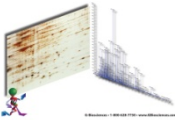
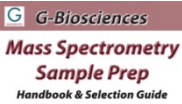
We recommend you wear gloves at all times and rinse them occasionally to reduce static build-up that attracts dust, hair and other interfering particles. Perform the entire process in a laminar flow hood, using tubes, tips and pipettes that were stored in the hood in a dust free environment. Avoid the use of detergents such as Triton and Tween (polymeric detergents) for cleaning flasks and glass plates used in electrophoresis.

PROTOCOL

1. Prepare fresh working SilverOUT™ reagents by mixing equal volumes SilverOUT™ I and II.
NOTE: For each protein band you will require ~50µl.
2. Add 50µl working SilverOUT™, ensuring the gel pieces are completely covered and vortex for 10 seconds. Incubate for 5-10 minutes or until the silver stain disappears from the gel band.
3. Remove working SilverOUT™ reagent and add 0.5ml ultrapure water, vortex and incubate for 5 minutes. Repeat the wash with ultrapure water until gel is clear.
4. Proceed with downstream applications.

RELATED PRODUCTS

Download our Mass Spectrometry Sample Preparation Handbook.



<http://info.gbiosciences.com/complete-mass-spectrometry-sample-preparation-handbook/>

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

Last saved: 11/20/2013 CMH



www.GBiosciences.com