



410PR

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

A Geno Technology, Inc. (USA) brand name

Protein-OUT™

For the Preparation of Plastic & Glassware
for Mass Spectrometry Analysis

(Cat. # 786-680)



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

INTRODUCTION

Protein-OUT™ is a unique washing solution that cleans plastic and glassware used in the proteolytic digestion and peptide isolation for mass spectrometry analysis. The rinsing of plastic and glassware with Protein-OUT™ will remove interfering chemicals and peptides that can leach from the tube walls and interfere or suppress the mass spectrometry signal. In addition, rinsing with Protein-OUT™ removes inadvertent contamination with dust and keratin. A simple rinse will prevent the need for repeat experiments, saving both time and money.

Protein-OUT™ is supplied in 50ml bottles and this is suitable for the rinsing of ~200 1.5-2ml micro centrifuge tubes.

ITEM(S) SUPPLIED (Cat. # 786-680)

Description	Size
Protein-OUT™	50ml

STORAGE CONDITION

Protein-OUT™ is shipped at ambient temperature. Upon arrival, store at room temperature.

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.

CAUTION

Protein-OUT™ is a volatile and corrosive product and should only be used as indicated below. Always wear gloves and safety glasses when using.

PROTOCOL

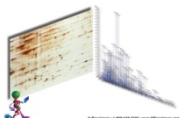
1. In a fume hood, open the Protein-OUT™ and pipette 250µl directly into the micro centrifuge tubes to be used.
NOTE: For larger sized containers use 250µl for every 2ml volume size. Ensure the container can be tightly sealed and then continue with the procedure.
2. Seal the tubes and vortex for 30 seconds to thoroughly rinse the tubes.
3. Remove the Protein-OUT™ from the tubes and discard in a disposal container containing >1L water. The diluted Protein-OUT™ can be disposed off down the sink.
4. Allow the tubes to dry completely before use.

RELATED PRODUCTS

Download our Mass Spectrometry Sample Preparation Handbook.



**Mass Spectrometry
Sample Prep**
Handbook & Selection Guide



<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: 10/4/2012 CMH



www.GBiosciences.com