



496PR

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

A Geno Technology, Inc. (USA) brand name

Immobilized Soybean Trypsin Inhibitor

For Removal of Trypsin, Chymotrypsin & Elastase

(Cat. # 786-843)



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

INTRODUCTION

G-Biosciences' Immobilized Soybean Trypsin Inhibitor is designed for the efficient removal of trypsin, chymotrypsin and elastase from protein digestions. The benefit of immobilizing the trypsin inhibitor is that there is no cross contamination of the purified samples with the soybean trypsin inhibitor, allowing for straightforward sample purifications.

An additional advantage is that the Immobilized Soybean Trypsin Inhibitor can be regenerated 10 times without significant loss in activity.

ITEM(S) SUPPLIED (Cat. # 786-843)

Description	Size
Immobilized Soybean Trypsin Inhibitor	2ml Resin

STORAGE CONDITION

It is shipped at ambient temperature. Upon arrival, store at -20°C.

PROPERTIES

- Binding Capacity: 9mg trypsin/ml resin
- Support: 4% Cross-linked Agarose

ADDITIONAL MATERIALS REQUIRED

- Disposable gravity flow or spin columns
www.gbiosciences.com/ResearchProducts/Protein-Research/Purification-and-Chromatography/Columns.aspx
- Binding Buffer (50mM Tris.HCl, 100mM NaCl, 10mM CaCl₂, pH 7.2)
- Elution Buffer (0.1M Acetic acid, 10mM CaCl₂)

PROTOCOL

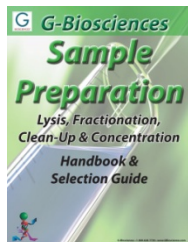
1. Prepare the protein sample by dialyzing or performing a buffer exchange with Binding Buffer.
NOTE: *We recommend our Tube-O-DIALYZER™ for dialysis of small samples and our SpinOUT™ columns for buffer exchange.*
2. Pack an appropriate volume of Immobilized Soybean Trypsin Inhibitor into a disposable column.
3. Equilibrate the resin with 5 column volumes (CV) of Binding Buffer.
4. Add the dialyzed protein sample to the column.
5. Slowly, apply Binding Buffer to the column and collect the flow-through, which contains the trypsin free protein sample.
6. The elution of the protein can be monitored by absorbance at 280nm. Apply Binding Buffer until the absorbances reach a baseline.
7. The bound trypsin can be eluted with the addition of Elution Buffer and monitoring the flow-through by absorbance at 280nm.

Column Regeneration

1. Ensure all bound trypsin has been eluted from the column by treating with Elution Buffer, until absorbances at 280nm reach a baseline.
2. Wash the column with 10 CV of Binding Buffer. The column can be regenerated 10 times without significant loss of activity.
3. For long term storage at -20°C, store in 50% glycerol in Binding Buffer supplemented with 0.05% sodium azide. Do not freeze unless stored in 50% glycerol.

RELATED PRODUCTS

Download our Sample Preparation Handbook



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

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

Last saved: 7/27/2012 CMH



www.GBiosciences.com