



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

A Geno Technology, Inc. (USA) brand name

Immobilized Trypsin

TPCK Treated Trypsin Immobilized On 4% Agarose

(Cat. #786-792)



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

INTRODUCTION

Immobilized Trypsin is TPCK treated Trypsin immobilized on 4% agarose that eliminates the contamination of protein digests by the trypsin. The immobilized trypsin is readily removed by separating the agarose from the digestion solution.

Trypsin is a serine endopeptidase that specifically cleaves peptide bonds on the carboxy side of *s*-aminoethyl cysteine, arginine and lysine residues and typically there is little or no cleavage at arginyl-proline and lysyl-proline bonds. The distribution of these residues in proteins allows trypsin digestion to produce peptides that are readily identified by mass spectrometry.

Native trypsin is prone to autolysis that results in pseudotrypsin, which exhibits a broader proteolytic specificity (a chymotrypsin like activity) and trypsin fragments that interfere with sequence analysis.

The Trypsin is TPCK treated to inactivate the interfering chymotrypsin activity and the resulting protein is affinity purified.

Immobilized Trypsin is supplied as a 50% slurry containing glycerol and sodium azide as a preservative.

ITEM(S) SUPPLIED (Cat. # 786-792)

Cat. #	Description	Size
786-792	Immobilized Trypsin	2ml resin

STORAGE CONDITIONS

Shipped at ambient temperature. Upon receipt store at 4°C, do NOT freeze.

IMPORTANT INFORMATION

Source: Bovine

Activity: ≥200 TAME units/ml resin (1 unit= 1μmole TAME (*p*-toluenesulfonyl-L-arginine methyl ester) hydrolyzed/min at pH8.2, 25°C)

Support: 4% Cross-linked Agarose

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.

ADDITIONAL COMPONENTS

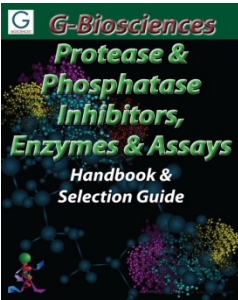
Digestion Buffer: 50mM ammonium bicarbonate, pH7.8

PROTOCOL

1. Prepare a 0.5ml 2mg/ml protein solution in 0.5ml Digestion Buffer.
2. Wash 100-250µl Immobilized Trypsin three times with 500µl Digestion Buffer.
3. Centrifuge the resin at 5,000rpm for 1-2 minutes.
4. Suspend the resin in 200µl Digestion Buffer and add the protein solution to the washed resin.
5. Incubate the reaction at 37°C in a shaking water bath for 2-18 hours.
6. Recover the digested protein solution by centrifuging at 5,000rpm for 1-2 minutes.

RELATED PRODUCTS

Download our Protease & Phosphatase Inhibitors, Enzymes & Assays Handbook.



<http://info.gbiosciences.com/protease-phosphatase-inhibitors-enzymes-assay-handbook>

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

Last saved: 11/4/2013 AT



www.GBiosciences.com