

Immobilized Soybean Trypsin Inhibitor

20235

1830.0

Number	Description
20235	Immobilized Soybean Trypsin Inhibitor , 2 ml settled gel, supplied as a 50% slurry in 50% glycerol and 0.05% sodium azide Binding Capacity: 9 mg of trypsin/ml of settled gel Support: 4% beaded agarose

Storage: Upon receipt store at -20°C. Product is shipped at ambient temperature.

Introduction

The Immobilized Soybean Trypsin Inhibitor is an effective tool for removing trypsin, chymotrypsin and elastase from protein digests. The soybean trypsin inhibitor is immobilized on beaded agarose enabling easy and efficient sample purifications. After purification, the support can be regenerated at least 10 times without significant loss in activity.

Procedure for Removing Trypsin from Samples

A. Materials Required

- Disposable gravity-flow or spin columns (for example, Disposable Polypropylene Columns, Product No 29922 or Handee™ Mini-Spin Columns, Product No. 69705)
- Binding Buffer: 50 mM Tris•HCl, 0.1 M NaCl, 10 mM CaCl₂; pH 7.2
- Protein sample dialyzed or desalted using the Binding Buffer
- Elution Buffer: 0.1 M acetic acid, 10 mM CaCl₂

B. Method

1. Pack a column with Immobilized Soybean Trypsin Inhibitor.
2. Equilibrate the column with Binding Buffer.
3. Add the dialyzed protein sample to the column.
4. Add Binding Buffer to the column and collect the flow-through, which contains the trypsin-free sample. Monitor sample collection by measuring its absorbance at 280 nm. Continue to add Binding Buffer until the absorbance approaches baseline.
5. To elute the bound trypsin, add Elution Buffer to the column and monitor flow-through by measuring its absorbance at 280 nm.
6. Regenerate the column by washing with Binding Buffer. The Immobilized Soybean Trypsin Inhibitor can be regenerated at least 10 times without significant loss in activity.
7. For storage, equilibrate column with 50% glycerol and 0.05% sodium azide and store at -20°C.

Note: The beaded agarose will become damaged if stored without 50% glycerol.

Related Pierce Products

78410	Halt™ Protease Inhibitor Cocktail Kit
78415	Halt™ Protease Inhibitor Cocktail, EDTA-Free
89834	Lysozyme, 25 g
89835	DNase I, 5,000 units
78115	Inclusion Body Solubilization Reagent, 100 ml
25200-25244	Precise™ Protein Gels (see catalog or web site for a complete listing)
24615	Imperial™ Protein Stain, 1 L, coomassie R-250 stain
24590	GelCode™ Blue Stain Reagent, 500 ml, coomassie G-250 stain
24582	E-Zinc® Reversible Stain Kit
26681	BlueRanger® Prestained Protein Molecular Weight Marker Mix
23227	BCA™ Protein Assay Kit
23236	Coomassie Plus – The Better Bradford™ Assay Kit

References

- Peterson, L. M., *et al.* (1976). Purification and crystallization of human carboxypeptidase A. *Biochemistry-USA*. **15(12)**: 2501-8.
- Reeck, G.R., *et al.* (1971). Isolation and characterization of carboxypeptidases A and B from activated pancreatic juice. *Biochemistry-USA*. **10(25)**:4690-8.
- Wu, C. *et al.* (1999). Proteolysis of native proteins. *J. Biol. Chem.* **274(2)**:1108-15.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current versions of product instructions are available at www.thermo.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2010 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.