

I. Introduction

This protocol is provided for **Capturem Trypsin** (Cat. No. 635722), single-use disposable mini spin columns containing membrane-immobilized trypsin for easy, rapid, and complete digestion of protein samples in less than 2 minutes. The columns are supplied together with an activation buffer.

II. Materials and Reagents

A. Components

- 20 Capturem Trypsin Columns
(mini spin columns containing a blue insert, supplied in 2-ml collection tubes)
- 5 ml Capturem 1X Activation Buffer

B. Additional Materials Required

1. Digestion Buffer

This product is compatible with commonly used ammonium bicarbonate and Tris-based buffers. Examples are shown below, but other buffers may also be used if necessary.

- 50–100 mM Tris, pH 8.0
- 10–50 mM ammonium bicarbonate, pH 8.0

2. Collection Tubes

Each sample will require 1 additional standard 2-ml collection tube, with or without a cap. These tubes should be used to collect peptides generated by digestion of protein samples which will be used for downstream analysis such as SDS-PAGE, HPLC, mass spectrometry, etc.

III. Sample Preparation

Some protein samples may require denaturation and/or reduction before enzymatic digestion. An example of a denaturation/reduction protocol is provided below. However, other protocols may be used if necessary.

1. Dissolve up to 80 µg of target protein or protein extract in your digestion buffer (see Section II.B.1) containing 6–8 M urea, in a reaction volume of 25–100 µl.
2. Add either DTT or TCEP to a concentration of 5 mM, vortex to mix, and incubate at 37°C for 30 minutes.
3. After denaturation, add a sufficient volume of digestion buffer (200–800 µl, as needed) to reduce the urea concentration to 1 M or below.

IV. Sample Digestion

Digest the protein samples prepared in Section III as follows:

1. Insert the Capturem Trypsin Column into the provided 2-ml collection tube.
2. Load 200 µl of Capturem 1X Activation Buffer onto the Capturem Trypsin Column to activate the column. Centrifuge at 500g for 1 min. Discard the flowthrough along with the collection tube and place the column in a new collection tube (supplied by the user—see Section II.B).
3. Load the protein sample from Section III, Step 3 onto the activated Capturem Trypsin Column. Centrifuge at 500g for 1 min. The eluted peptides are ready for downstream analysis.

NOTE: Capturem Trypsin digestion may generate longer and more unique peptides than in-solution trypsin digestion, due to an increase in missed cleavage sites. Identification of these unique peptides may require adjustments to the minimum/maximum peptide mass and the maximum number of allowed missed cleavages in bioinformatics data analysis, to obtain complete sequence coverage of the target protein.

| Contact Us | |
|--|--|
| Customer Service/Ordering | Technical Support |
| tel: 800.662.2566 (toll-free) | tel: 800.662.2566 (toll-free) |
| fax: 800.424.1350 (toll-free) | fax: 800.424.1350 (toll-free) |
| web: takarabio.com | web: takarabio.com |
| e-mail: ordersUS@takarabio.com | e-mail: techUS@takarabio.com |

Notice to Purchaser

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Takara Bio USA, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

©2017 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

This document has been reviewed and approved by the Quality Department.