



239PR

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

A Geno Technology, Inc. (USA) brand name

Trypsin Digestion Mix

Provides optimal buffered conditions for in gel trypsin digestion of proteins

(Cat. #786-242)



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

INTRODUCTION 3

ITEM(S) SUPPLIED (Cat. # 786-242) 3

STORAGE CONDITION 3

Important Information 3

Protocol 4

 A. Excise Protein Spots/Bands 4

 B. In-Gel Reduction, Alkylation & Destaining of Proteins 4

 Silver Destain, Denaturation and Alkylation 4

 C. In Gel Trypsin Digestion 5

 Preparation Before Use 5

 D. Peptide Extraction 6

 For MALDI peptide mass mapping..... 6

 For LC MS/MS 6

Related Products 7

INTRODUCTION

The following protocol is a generic protocol for the in gel digestion of proteins in a silver stained gel using our Trypsin Digestion Mix. For optimal in gel digestion, we recommend our InGel™ Silver kit (Cat. # 786-241).

The first section of the protocol covers the in-gel digestion of proteins by trypsin. Trypsin is a serine endopeptidase that specifically cleaves peptide bonds on the carboxy side of s-aminoethyl cysteine, arginine and lysine residues. Typically there is little or no cleavage at arginyl-proline and lysyl-proline bonds.

The second part of the protocol focus on the extraction of the digested proteins from the gel pieces for analysis by mass spectrometry (MALDI and LC MS/MS)

ITEM(S) SUPPLIED (Cat. # 786-242)

Description	Size
Trypsin Digestion Mix	50ml

STORAGE CONDITION

Shipped at ambient temperature. Store at RT.

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.

Reduction & Alkylation will minimize artifactual peaks caused by disulfide bridges and side chain modifications and improve detection of peptides with cysteines. Alkylation by iodoacetamide will increase the mass of peptides by 57.02/cysteine present.

PROTOCOL

A. Excise Protein Spots/Bands

Processing of protein bands/spots. Following electrophoresis the proteins need to be fixed in the gel matrix. If a fixing step is not included with your silver staining technique we recommend fixing in 5% acetic acid in 1:1 ultrapure water: methanol. For silver staining avoid the use of cross-linking reagents (i.e. glutaraldehyde) or strong oxidizers (i.e. chromates or permanganates). We recommend FOCUS™ FASTsilver™ (Cat. # 786-240), a mass spectrometry compatible silver stain.

1. Rinse the entire gel in ultrapure water for 1-2 hours before processing.
2. Excise protein spots or bands with a clean scalpel and cut bands to 1-2mm cubes.

***NOTE:** Pieces smaller than 1mm² may clog pipette tips in further processing.*

3. Transfer to 0.5ml clean centrifuge tubes and briefly spin down in a benchtop centrifuge.

***NOTE:** Tubes can be cleaned with Protein-OUT™ (Cat. # 786-680), a unique solution to remove proteins and other mass spectrometry interfering agents.*

B. IN-GEL REDUCTION, ALKYLATION & DESTAINING OF PROTEINS

Silver Destain, Denaturation and Alkylation

1. Destain the gel piece with ~50µl silver destain, ensuring the gel pieces are completely covered and vortex for 10 seconds. Incubate for 5-10 minutes or until the silver stain disappears from the gel band. We recommend SilverOUT™ (Cat. # 786-244)
2. Remove the destain reagent and add 0.5ml ultrapure water, vortex and incubate for 5 minutes. Repeat the wash with ultrapure water until gel is clear.
3. Add 500µl acetonitrile and incubate at room temperature for 10 minutes, or until the gel pieces become opaque and shrink. Briefly centrifuge to pellet the gel pieces and remove all the liquid.
4. Add 20-50µl 10mM DTT solution, ensuring the gel pieces are completely covered. Incubate at 60°C for 30 minutes.
5. Allow the tubes to cool to room temperature and then repeat step 4.
6. Add 20-50µl 50mM iodoacetamide solution, ensuring the gel pieces are completely covered. Incubate at room temperature for 20 minutes in the dark.

7. Add 500µl acetonitrile and vortex and incubate for 5 minutes. Discard the acetonitrile.
8. Repeat the acetonitrile wash (step 8) until the gel pieces are opaque white and completely dehydrated. This normally requires 2-3 washes.
9. Dry the gel pieces in a vacuum centrifuge and store at -20°C until use.

C. IN GEL TRYPSIN DIGESTION

Preparation Before Use

Trypsin Rehydration Solution: To 1.2ml ultrapure water, add 150µl Trypsin Digestion Mix and 150µl acetonitrile and vortex.

Trypsin Working Solution: Add 1.5ml Trypsin Rehydration Solution to 20µg mass spectrometry grade trypsin. We recommend our Trypsin, Mass Spectrometry Grade (Cat. # 786-245). Incubate for 15-30 minutes on ice and then gently pipette up and down to fully rehydrate. Do not vortex as this will lead to a loss of activity. The Trypsin should be solubilized immediately before use and unused Trypsin discarded.

***NOTE:** If smaller volumes of trypsin are required, resuspend in an appropriate volume of Trypsin Suspension Buffer, incubate for 15-30 minutes on ice and then gently pipette up and down to fully rehydrate. Aliquot into 10µl aliquots and store at -20°C. To use, add 150µl Trypsin Rehydration Solution to each 10µl fraction.*

1. Add ~50µl Trypsin Working Solution to the gel pieces (enough to cover the gel pieces)
2. Incubate at 4°C for 30 minutes.
3. If all the Trypsin Working Solution has been absorbed then add more Trypsin Working Solution, ensuring the gel pieces remain covered.
4. Incubate for a further 90 minutes at 4°C and then add 10-20µl Trypsin Digestion Mix to cover the gel pieces.
5. Incubate the digestion tube at 37°C for overnight for maximal peptide recovery.

D. PEPTIDE EXTRACTION

For MALDI peptide mass mapping

1. Cool the tubes to room temperature, centrifuge briefly in a benchtop centrifuge and remove a 1.5µl aliquot for MALDI peptide mapping.

For LC MS/MS

1. Briefly, centrifuge the digestion and add 40µl *Pep-Extract*[™] (Cat. # 786-243) and vortex. Incubate at 37°C for 15-30 minutes with periodic vortexing.
2. Centrifuge the tube briefly and collect the extract for analysis using a fine tip pipette to prevent removing the fine-gel particles that may clog analysis equipment. If samples require guanidation continue to the next section.
3. Dry the extra sample in a vacuum centrifuge and store at -20°C until use. Dried extracts can be stored for a few months.
4. To reconstitute for further LC MS/MS, add 10µl 0.1% trifluoroacetic acid and vortex. Remove the required aliquot for analysis and then dry the remaining sample in a vacuum centrifuge and store at -20°C.
5. The samples can now be guanidated.

RELATED PRODUCTS

Download our Mass Spectrometry Handbook.



Mass Spectrometry

Sample Prep

Handbook & Selection Guide



<http://info2.gbiosciences.com/complete-mass-spectrometry-sample-preparation-handbook>

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

Last saved: 6/10/2015 CMH



www.GBiosciences.com