## **INSTRUCTIONS**



# Pierce LysN Protease, MS Grade

90300 90301

Number Description

90300 Pierce LysN Protease, MS Grade, 20μg
 90301 Pierce LysN Protease, MS Grade, 5 × 20μg

**Storage:** Upon receipt, store at -20°C. Product is shipped on ice packs.

#### **Table of Contents**

Table o	of Contents	. 1		
Introdu	Introduction			
Important Product Information				
Material Preparation				
Procedure for In-solution Protein Digestion				
	Additional Materials Required			
	Reduction and Alkylation			
	Digestion			
Troubleshooting				
Related Thermo Scientific Products				
	General References			

#### Introduction

Thermo Scientific<sup>TM</sup> Pierce<sup>TM</sup> LysN Protease is a mass spectrometry (MS)-grade metalloprotease that cleaves at the aminoside of lysine residues. LysN is a zinc metalloprotease derived from *Grifola frondosa* that is resistant to moderate denaturing conditions making it ideal for digestion of samples for "shotgun" proteomics. Peptides generated by LysN are longer than those generated by trypsin and have more prevalent charged amino-terminal peptide fragments which are useful for *de novo* sequencing with collision induced dissociation (CID). The purified native Pierce LysN Protease has improved stability, specific activity and cleavage selectivity.

### **Important Product Information**

- Maximal LysN activity occurs at pH 7-9. Common digestion buffers include 50mM ammonium bicarbonate, pH 8;
   50mM Tris, pH 8; and 50mM TEAB, pH 8.5.
- LysN remains active under moderate denaturing conditions including 0.1% SDS, 6M urea or heat up to 70°C.
- EDTA or other metal chelators interfere with LysN activity. To recover activity loss due to the presence of metal chelators, add a zinc salt such as ZnSO<sub>4</sub> at a concentration greater than the metal chelator to the digestion buffer.
- Reconstituted stock solutions of LysN in water are stable at -80°C for 2 years or -20°C for one year without significant loss in activity. Minimize the number of freeze/thaw cycles by aliquoting stock solutions of the enzyme.
- Reduction and alkylation of cysteine residues using dithiothreitol (DTT) and iodoacetamide (IAA), respectively, will
  cleave disulfide bonds and prevent disulfide bond reformation. These steps improve digestion of cysteine-containing
  proteins and detection of cysteine-containing peptides. Alkylation with IAA increases the mass of a peptide by 57.02Da
  for each cysteine present.



#### **Material Preparation**

Enzyme Reconstitution Reconstitute Pierce LysN Protease using LC/MS grade water to 0.5 mg/mL (i.e., add  $40 \mu L$ 

LC/MS grade water to 20µg of lyophilized LysN). Aliquot reconstituted enzyme in single-use

volumes and store at -20°C or -80°C.

## **Procedure for In-solution Protein Digestion**

Note: The following protocol is an example application for this product. Specific applications will require optimization.

#### A. Additional Materials Required

- 1M Tris, pH 8 (e.g., Fisher Scientific Product No. BP1758-100)
- Urea, sequanal grade (e.g., Thermo Scientific Product No. 29700)
- Ammonium bicarbonate (e.g., Acros Product No. 370930250)
- DTT (e.g., Thermo Scientific Product No. 20290 or 20291)
- IAA (e.g., Thermo Scientific Product No. 90034)
- LC/MS grade water (e.g., Thermo Scientific Product No. 51140)
- Optional: 0.5M TCEP (e.g., Thermo Scientific Product No. 77720)
- Optional: SDS (e.g., Fisher Scientific Product No. BP1311-1)
- Optional: Thermo Scientific<sup>TM</sup> Pierce<sup>TM</sup> C18 Spin Columns (Product No. 89870)

#### B. Reduction and Alkylation

 Dissolve protein in 50mM ammonium bicarbonate or 50mM Tris, pH 7.5-8 with a denaturing reagent such as 8M urea or 0.1% SDS.

**Note:** Sodium deoxycholate is not recommended as a denaturing reagent because it reduces LysN activity.

- 2. Prepare a fresh solution of 500mM DTT by dissolving 7.7mg of DTT in 100μL of ultrapure water.
- 3. Add 500mM DTT solution to the protein sample to a final concentration of 20mM (1:25 dilution) and mix briefly.
- 4. Incubate at 37°C for 45 minutes or 95°C for 10 minutes.
- 5. Prepare a fresh solution of 1M IAA by dissolving 93mg of IAA in 500μL of ultrapure water.

Note: Protect IAA stock solutions from light.

- 6. Add 1M IAA solution to the reduced protein sample to a final concentration of 40mM (1:25 dilution) and mix briefly.
- 7. Incubate the reaction mixture at room temperature for 15 minutes protected from light.
- 8. Quench the alkylation reaction by adding 500mM DTT solution to a final concentration of 10mM (1:50 dilution).

#### C. Digestion

 $1. \quad \text{Add reconstituted LysN solution to the sample to a final protease-to-protein ratio of 1:75 to 1:100 (w/w).} \\$ 

**Note:** Protein samples dissolved in 8M urea must be diluted to  $\leq$  6M urea before digestion. For SDS-containing samples, dilution is not necessary.

2. Incubate the sample at 50°C for 2 hours or 37°C for 4 hours.

**Note:** Digestion can also be performed at 70°C for 45 minutes; however, more missed cleavages will occur.

3. Store samples at -20°C to stop digestion reactions. Immediately before MS analysis, clean-up samples with C18 spin columns (Pierce C18 Spin Columns, Product No 89870) ensuring to acidify the sample with formic acid for proper C18 binding.



# **Troubleshooting**

Problem	Possible Cause	Solution
No or incomplete digestion	Incorrect pH or buffer conditions	Check buffer pH
	Reduced enzymatic activity	Once reconstituted, store enzyme at -80°C for 1 year. Do not let enzyme sit at room temperature
		Ensure no metal chelator is present or add a Zn <sup>2+</sup> salt
	Too little enzyme used	Increase amount of LysN for digestion
	Too high incubation temperature used	Decrease incubation temperature
Low sequence coverage	Too few, too many or unevenly distributed protease digestion sites	Separately use multiple proteases to digest the sample and combine results (e.g., multi-consensus reports in Thermo Scientific <sup>TM</sup> Proteome Discoverer Software)
Precipitation after alkylation	Too much reduction/alkylation buffer for quantity of protein being digested	Quench alkylation reaction using 10mM DTT
Over-alkylation	Alkylation was allowed to proceed for too long	Alkylate at room temperature for 30 minutes and quench reaction with 10mM DTT
Incomplete alkylation or incomplete recovery of alkylated peptides	Used old or inactive iodoacetamide solution	Prepare iodoacetamide solution immediately before use, and protect it from light
Too much background noise during LC-MS	Buffer, salt or urea interference	Clean-up sample before analysis with reversed-phase tips or spin cartridges (e.g., Pierce C18 Spin Columns)

## **Related Thermo Scientific Products**

90057	Pierce Trypsin Protease, MS Grade, $5 \times 20 \mu g/vial$
90058	Pierce Trypsin Protease, MS Grade, $5 \times 100 \mu g/vial$
90059	Pierce Trypsin Protease, MS Grade, 1mg/vial
84849	Pierce Mass Spec Sample Prep Kit for Cultured Cells
20233	Immobilized TPCK Trypsin, 50mg
90051	Lys-C Endoproteinase, MS Grade, $20\mu g$
90054	Glu-C Endoproteinase, MS Grade, $5 \times 10 \mu g$
90053	Asp-N Endoproteinase, MS Grade, 2μg
90056	Chymotrypsin Endoproteinase, TLCK treated, MS Grade, $4\times25\mu g$
89895	In-Solution Tryptic Digestion and Guanidination Kit
89870	Pierce C18 Spin Columns, 25/pkg
28904	Trifluoroacetic Acid, Sequanal Grade, $10 \times 1 \text{mL}$



**28905** Formic Acid,  $10 \times 1$ mL

88300 Pierce Fe-NTA Phosphopeptide Enrichment Kit

88301 Pierce TiO<sub>2</sub> Phosphopeptide Enrichment and Clean-up Kit

**88302** Pierce Graphite Spin Columns, 30 columns

#### **General References**

Taouatas, N., et al. (2010). Evaluation of metalloendopeptidase Lys-N protease performance under different sample handling conditions. J Proteome Res. 9(8):4282-8.

Gershon, P. (2014). Cleaved and missed sites for trypsin, Lys-C, and Lys-N can be predicted with high confidence on the basis of sequence content. *J Proteome Res.* **13(2)**:702-9.

Nonaka, T. J., et al. (1998). Kinetic characterization of lysine-specific metalloendopeptidases from *Grifola frondosa* and *Pleurotus ostreatus* fruiting bodies. *Biochemistry* 124(1):157-62.

Products are 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"). 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 Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

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 REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR 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.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at <a href="https://www.thermoscientific.com/pierce">www.thermoscientific.com/pierce</a>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2014 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.