

## Certificate of Analysis

### Elastase:

**Part No.** V189A  
**Size** 5mg

**Description:** Elastase is a serine protease that preferentially cleaves at the C-terminus of alanine, valine, serine, glycine, leucine or isoleucine (1–4). Elastase has a unique capability of digesting elastin (5). This enzyme can be used alone or in combination with other proteases for protein analysis by mass spectrometry and other applications.

**Biological Source:** Porcine pancreas.

**Molecular Weight:** 25.9kDa (6).

**Form:** Lyophilized.

**Storage Conditions:** See the Product Information Label for storage conditions and expiration date.

**Optimal pH:** 9.0.

**Activators:** Elastase is activated by sodium carbonate, sodium sulfate and Tris (7).

#### Inhibitors:

- Irreversible:**  $\beta$ -casomorphin-7 (BCM7) (5); pH 3–4 (8); diisopropyl-phosphofluoridate and alkyl isocyanates (9); peptide chloromethyl ketone (10,11).
- Competitive:** Derivatives of dipeptides and alanine, valine, leucine and isoleucine (12).
- Selective:** Soybean trypsin inhibitor and kallikrein inhibitor suppress proteolytic but not elastolytic activity (13).

**Usage Note:** Resuspend Elastase in double-distilled water to a final concentration of 1mg/ml. Store reconstituted Elastase at 4°C for up to 2 weeks.

## Quality Control Assays

This lot passes the following Quality Control specifications:

**Activity:** Digestion reactions using glucagon as a substrate at either a 1:20 or 1:100 protease:substrate ratio show no detectable intact substrate remaining by reverse-phase HPLC analysis after 30 minutes of digestion at 37°C.

### Usage Information on Back

Signed by:



R. Wheeler, Quality Assurance

Part# 9PIV189

Revised 8/16



AF9PIV189 0816V189



## Promega

### Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	<a href="http://www.promega.com">www.promega.com</a>

#### PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 2011–2013, 2016 Promega Corporation. All Rights Reserved.

ProteaseMAX is a trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIV189  
Printed in USA. Revised 8/16.

## 1. In-Solution Digestion Protocol

1. Resuspend the protein in reaction buffer.  
**Note:** Tris is an activator of Elastase and must be included in the reaction buffer.
2. Resuspend the Elastase in double-distilled water.
3. Transfer the protein solution to a microcentrifuge tube.
4. Add Elastase to protein solution; mix. We recommended using enzyme:protein ratios of 1:20 to 1:100.
5. Incubate 2–18 hours at 37°C.
6. Stop the reaction by adding 10% formic acid or TFA to a final concentration of 0.5% or by heating at 95°C for 10 minutes.

## 2. Composition of Buffers and Solutions

### reaction buffer

50mM Tris-HCl (pH 8.5–9.5)

## 3. References

1. Narayanan, S. *et al.* (1969) *Biochemistry J.* **114**, 11–7.
2. Gertler, A. *et al.* (1977) *Biochemistry* **16**, 2709.
3. Del Mar, E.G. (1980) *Biochemistry* **19**, 468–72.
4. Powers, J.C. *et al.* (1977) *Biochimica et Biophysica Acta* **485**, 156–66.
5. Wright, P.A. *et al.* (2001) *Eur J. Biochem* **268**, 2969–74.
6. *Handbook of Proteolytic Enzymes* (1998) 42–46.
7. Shotton, D.M. (1970) *Methods Enzymol.* **19**, 113–40.
8. Wasi, S. *et al.* (1968) *Biochem. J.* **106**, 926–7.
9. Brown, W. *et al.* (1973) *Biochemistry* **12**, 835–40.
10. Thompson, R.C. *et al.* (1973) *Biochemistry* **12**, 44–7.
11. Powers, J. *et al.* (1972) *J. Am. Chem. Soc.* **94**, 6544–5.
12. Dzialoszynski, L. *et al.* (1973) *Biochim Biophys Acta.* **302**, 406–10.
13. Walford, R.L. *et al.* (1962) *Arch. Biochem Biophys.* **98**, 191–6.

## 4. Related Products

Product	Size	Conc.	Cat. #
Asp-N, Sequencing Grade	2µg		V1621
Arg-C, Sequencing Grade	10µg		V1881
Chymotrypsin, Sequencing Grade	25µg		V1061
	100µg (4 × 25µg)		V1062
Endo H	10,000u	500u/µl	V4871
	50,000u	500u/µl	V4875
Endoproteinase Lys-C, Sequencing Grade	5µg		V1071
Fetuin	500µg	10mg/ml	V4961
Glu-C, Sequencing Grade	50µg (5 × 10µg)		V1651
Immobilized Trypsin	2ml		V9012
	4ml (2 × 2ml)		V9013
Pepsin	250mg		V1959
PNGase F	500u	10u/µl	V4831
ProteaseMAX™ Surfactant, Trypsin Enhancer	1mg		V2071
	5 × 1mg		V2072
Protein Deglycosylation Mix	20 reactions		V4931
rLys-C, Mass Spec Grade	15µg		V1671
Sequencing Grade Modified Trypsin	100µg (5 × 20µg)		V5111
Sequencing Grade Modified Trypsin, Frozen	100µg (5 × 20µg)		V5113
Thermolysin	25mg		V4001
Trypsin Gold, Mass Spectrometry Grade	100µg		V5280
Trypsin/Lys-C Mix, Mass Spec Grade	20µg		V5071
	100µg		V5072
	100µg (5 × 20µg)		V5073