

Certificate of Analysis

PNGase F

Part No.	Size
V483A	500u

Description: N-Glycosidase F (PNGase F) catalyzes the cleavage of N-linked oligosaccharides. PNGase F is a recombinant glycosidase cloned from *Elizabethkingia miricola*.

Biological Source: *E. coli*.

Concentration: 10,000u/ml.

Molecular Weight: PNGase F has a molecular weight of approximately 36kDa.

Physical Form: PNGase F is supplied as a liquid in 20mM Tris-HCl (pH 7.5 at 25°C), 50mM NaCl and 5mM EDTA at a concentration of 10,000u/ml.

Storage Conditions: Store at +2° to +10°C.

Unit Definition: One unit of PNGase F will catalyze the deglycosylation of 1 nanomole of denatured Ribonuclease B (RNase B) in one minute at 37°C. One Promega unit is equal to 1 IUB milliunit.

Part# 9PIV483

Revised 1/18



AF9PIV483 0118V483

Quality Control Assays

This lot passes the following Quality Control specifications:

Activity Assay: Denatured RNase B (20µg) is incubated with PNGase F for 10 minutes at 37°C and then analyzed by SDS-PAGE. Fully glycosylated RNase B migrates at approximately 17kDa. Deglycosylation is assessed by the presence of deglycosylated RNase B with an apparent molecular weight of 13.7kDa.

Purity: ≥95% as determined by SDS-PAGE analysis.

Usage Information on Back



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	www.promega.com

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.

© 2012–2013, 2016, 2018 Promega Corporation. All Rights Reserved.

ProteaseMAX is a trademark of Promega Corporation. ZipTip is a registered trademark of Millipore Corporation. Triton is a registered trademark of Union Carbide Chemicals & Plastics Technology 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# 9PIV483
Printed in USA. Revised 1/18

Signed by:

R. Wheeler, Quality Assurance

1. Protein Deglycosylation Using Recombinant PNGase F

Note: The following protocols are intended as a general guide for protein deglycosylation. Activity against different glycoprotein substrates is highly dependent on reaction conditions and should be determined empirically.

A. Protocol 1: Protein Deglycosylation Using Denaturing Conditions for SDS-PAGE

Materials to Be Supplied By the User

- 5% SDS
- 1M DTT
- 0.5M sodium phosphate buffer (pH 7.5)
- 10% NP-40

1. Add up to 50µg of the target glycoprotein in water (or a compatible buffer at a low ionic strength) to a final volume of 12µl.
2. Add 1µl of 5% SDS.
3. Add 1µl of 1M DTT.
4. Denature sample by heating at 95°C for 5 minutes.
5. Cool sample for 5 minutes at room temperature.
6. Add 2µl of 0.5M sodium phosphate buffer (pH 7.5).
Note: Other buffers can be used if they are within the acceptable pH range for PNGase F (pH 6–10).
7. Add 2µl of 10% NP-40.
Note: Triton® X-100 may be substituted for NP-40.
8. Add 2µl of recombinant PNGase F.
9. Incubate at 37°C for 1–3 hours.

Notes:

Deglycosylation of glycoproteins may be visualized by gel-shift on SDS-PAGE, with the deglycosylated product running faster than the glycosylated substrate.

Samples produced by SDS-PAGE are compatible with mass spectrometry analysis following standard protocols. (Refer to the *Trypsin Gold, Mass Spectrometry Grade, Technical Bulletin*, #TB309).

B. Protocol 2: Protein Deglycosylation Using Non-Denaturing Conditions for Mass Spectrometry

Materials to Be Supplied By the User

- 50mM ammonium bicarbonate buffer (pH 7.8)
1. Add up to 20µg of glycoprotein in 50mM ammonium bicarbonate (pH 7.8) to a final volume of 18µl.
 2. Add 2µl of recombinant PNGase F.
 3. Incubate at 37°C for 2–18 hours.

Notes:

Most substrates are deglycosylated more effectively when denatured. Deglycosylation under nondenaturing conditions may require increasing both the amount of PNGase F used and the incubation time. Deglycosylation of some substrates may be enhanced by the addition of up to 0.1% ProteaseMAX™ Surfactant.

Samples are compatible with downstream mass spectrometry analysis using either gel- or solution-based digestion protocols. To desalt the sample prior to MS analysis, see the ZipTip® protocol provided in the *Trypsin Gold, Mass Spectrometry Grade, Technical Bulletin* #TB309.

2. References

1. *Trypsin Gold, Mass Spectrometry Grade, Technical Bulletin* #TB309, Promega Corporation.
2. Mann, A.C., Self, C.H., and Turner, G.A (1994) A general method for the complete deglycosylation of a wide variety of serum glycoproteins using peptide-N-glycosidase-F. *Glycoconjugate Journal* **1**, 253–61.

3. 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
Elastase	5mg		V1891
Endo H	10,000u	500u/µl	V4871
	50,000u	500u/µl	V4875
Fetuin	500µg	10mg/ml	V4961
Glu-C, Sequencing Grade	50µg (5 × 10µg)		V1651
Immobilized Trypsin	2ml		V9012
	4ml (2 × 2ml)		V9013
Pepsin	250mg		V1959
ProteaseMAX™ Surfactant, Trypsin Enhancer	1mg		V2071
	5 × 1mg		V2072
rLys-C, Mass Spec	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