PNGase F Glycan Cleavage Kit

Catalog Number A39245

Pub. No. MAN0018620 Rev. B

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

The PNGase F Glycan Cleavage Kit includes all components necessary to perform the enzymatic removal of almost all N-linked oligosaccharides from glycoproteins. The kit includes recombinant Peptide N-Glycosidase F (PNGase F) enzyme, which cleaves N-glycan chains at the innermost GlcNAc and asparagine residues of high mannose, hybrid, and complex oligosaccharides, and a 10X reaction buffer.

Contents and storage

Content	Amount	Storage ^[1]
Recombinant, PNGase F (500,000 units ^[2])	60 µL	–20°C
PNGase F 10X Buffer	1 mL	

^[1] Stable for up to 24 months.

 $^{2]}$ One unit is defined as the amount of enzyme required to remove > 95% of the carbohydrate from 10 µg of denatured RNase B in 1 hour at 37°C in a total reaction volume of 10 µL.

General guidelines

- Do not use SDS in the reaction mixture because SDS inhibits PNGase F activity. If detergent is needed, use NP-40.
- N-linked glycans containing core α1-3 Fucose are not cleaved by PNGase F.
- For deglycosylation of native glycoprotiens, increased incubation time and increased amount of enzyme may be needed.
- Optimize reaction conditions for each substrate.
- If a larger amount of glycoprotein is used, scale up reaction volumes accordingly.

Digest glycoprotein with PNGase F (nondenaturing conditions)

Note: The following instructions are for one sample. The Master Mix volume can be scaled up proportionally for multiple samples.

- 1. Prepare a Master Mix by combining 4 μ L of PNGase F 10X Buffer with 0.5 μ L of PNGase F enzyme.
- 2. Add 25 μ g of glycoprotein sample to a 1.5-mL tube, then bring the reaction volume up to 35.5 μ L with HPLC-grade water.
- 3. Add 4.5 μL of the prepared Master Mix, then mix gently.
- 4. Incubate the reaction at 50°C for 1 hour, then centrifuge briefly.
- 5. Analyze by method of choice.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/ home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.





Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, California 92008 USA For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN0018620 B

Revision	Date	Description
В	7 March 2025	Updated Recombinant, PNGase F amount in Contents and storage.
A.0	14 April 2019	New document.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019-2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.