

Product Information

Endoglycosidase F1
Catalog Number **E9762**

Endoglycosidase F2
Catalog Number **E0639**

Endoglycosidase F3
Catalog Number **E2264**

Storage Temperature 2–8 °C

TECHNICAL BULLETIN

Product Description

Endoglycosidases F1, F2, and F3 are intended for the deglycosylation of N-linked oligosaccharides from glycoproteins under native conditions. Particular residues, due to their location in the native protein structure, are often resistant to the traditional deglycosylation methods using PNGase F and cannot be removed unless the protein is denatured.

Endoglycosidases F1, F2, and F3 are less sensitive to protein conformation than PNGase F and are more suitable for deglycosylation of native proteins. The linkage specificities of Endoglycosidases F1, F2, and F3 suggest a general strategy for deglycosylation of proteins that may remove all classes of N-linked oligosaccharides without denaturing the protein (see Figure 1). All complex oligosaccharides can be reduced to the trimannosylchitobiose core by treatment of the glycoproteins with neuraminidase, β -galactosidase, and N-acetylglucosaminidase, available as part of the EDEGLY Kit (see Figure 2). Fucosidases may be required in some situations. The remaining trimannosylchitobiose core structures can be removed with Endoglycosidase F3. Biantennary and triantennary structures can be immediately removed by Endoglycosidases F2 and F3, respectively. Oligomannose and hybrid structures can be removed by Endoglycosidase F1.

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

These products ship on wet ice and storage at 2–8 °C is recommended. The products may be used for at least 1 year when stored as indicated. Several days exposure to ambient temperatures will not reduce the activity of the enzymes.

Endoglycosidase F1
from *Elizabethkingia miricola*
recombinant, expressed in *E. coli*
Catalog Number **E9762**

EC 3.2.1.96

Synonyms: Endo- β -N-acetylglucosaminidase F1,
Endo F1

Elizabethkingia miricola was formerly known as *Elizabethkingia*, *Chryseobacterium*, or *Flavobacterium meningosepticum*.

The type of carbohydrate structure, oligomannose, hybrid, or complex, and the state of $\alpha(1\rightarrow6)$ core fucosylation have profound effects on the cleavage by the endoglycosidases. Endoglycosidase F1 cleaves asparagine-linked or free oligomannose and hybrid, but not complex, oligosaccharides. Core fucosylation of hybrid structures reduces the rate of cleavage by Endo F1 more than 50-fold. The core fucosylated hybrids are usually only found on glycoproteins from swainsonine-treated cells. Endo F1 will cleave sulfated high-mannose oligosaccharides; whereas, Endo H will not. Endo F1 cleaves between the two N-acetylglucosamine residues in the diacetylchitobiose core of the oligosaccharide, generating a truncated sugar molecule with one N-acetylglucosamine residue remaining on the asparagine. In contrast, PNGase F removes the oligosaccharide intact (see Figure 3).

The enzyme has a molecular mass of ~32 kDa.

The enzyme is provided as a solution in 20 mM Tris-HCl, pH 7.5.

Deglycosylation Procedure

1. Add up to 200 µg of glycoprotein to a microcentrifuge tube. Adjust the final volume to 37.5 µl with deionized water.
2. Add 10 µl of 5× Reaction Buffer (Catalog Number R9025)
3. Add 2.0 µl of Endo F1 solution to the reaction. Incubate 1 hour at 37 °C.
4. Monitor cleavage by SDS-PAGE.

Endoglycosidase F2
from *Elizabethkingia miricola*
recombinant, expressed in *E. coli*
 Catalog Number **E0639**

EC 3.2.1.96

Synonyms: Endo-β-N-acetylglucosaminidase F2, Endo F2

Elizabethkingia miricola was formerly known as *Elizabethkingia*, *Chryseobacterium*, or *Flavobacterium meningosepticum*.

Endo F2 and Endo F3 are unique in their ability to cleave complex structures. Endo F2 cleaves asparagine-linked or free oligomannose, and biantennary complex oligosaccharides. Oligomannose structures are cleaved at a 20-fold reduced rate. Fucosylation has little effect on Endo F2 cleavage of biantennary structures. It will not cleave hybrid structures. It cleaves between the two N-acetylglucosamine residues in the diacetylchitobiose core of the oligosaccharide, generating a truncated sugar molecule with one N-acetylglucosamine residue remaining on the asparagine. Endo F2 is less sensitive to protein conformation than PNGase F and is more suitable for deglycosylation of native proteins (see Figure 4).

The enzyme has a molecular mass of ~32 kDa.

The enzyme is provided as a solution in 10 mM sodium acetate and 25 mM sodium chloride, pH 4.5.

Deglycosylation Procedure

1. Add up to 200 µg of glycoprotein to a microcentrifuge tube. Adjust the final volume to 37.5 µl with deionized water.
2. Add 10 µl of 5× Reaction Buffer (Catalog Number R9150)
3. Add 2.0 µl of Endo F2 solution to the reaction. Incubate 1 hour at 37 °C.
4. Monitor cleavage by SDS-PAGE.

Endoglycosidase F3
from *Elizabethkingia miricola*
recombinant, expressed in *E. coli*
 Catalog Number **E2264**

EC 3.2.1.96

Synonyms: Endo-β-N-acetylglucosaminidase F3, Endo F3

Elizabethkingia miricola was formerly known as *Elizabethkingia*, *Chryseobacterium*, or *Flavobacterium meningosepticum*.

Endo F3 is unique in that its cleavage is sensitive to the state of peptide linkage of the oligosaccharide, as well as the state of core fucosylation. Endoglycosidase F3 cleaves asparagine-linked biantennary and triantennary complex oligosaccharides. It will cleave non-fucosylated biantennary and triantennary structures at a slow rate, but only if peptide-linked. Core fucosylated biantennary structures are efficient substrates for Endo F3, even as free oligosaccharides. Endo F3 will also cleave fucosylated trimannosyl core structures on free and protein-linked oligosaccharides. Core fucosylation of biantennary structures increases activity up to 400-fold. There is no activity on oligomannose and hybrid molecules. It will only cleave non-fucosylated trimannosyl core structures when attached to a protein, not as free oligosaccharides. This activity was previously attributed only to Endoglycosidase D from *Diplococcus (Streptococcus) pneumonia* (see Figure 4). The effect of fucosylation on triantennary structures has not yet been investigated. Endo F3 cleaves between the two N-acetylglucosamine residues in the diacetylchitobiose core of the oligosaccharide, generating a truncated sugar molecule with one N-acetylglucosamine residue remaining on the asparagine. In contrast, PNGase F removes the oligosaccharide intact. When used in conjunction with the EDEGLY Kit, all complex oligosaccharides can be reduced to the trimannosylchitobiose core with subsequent removal with Endo F3.

Endo F3 is less sensitive to protein conformation than PNGase F and is, therefore, more suitable for deglycosylation of native proteins.

The use of detergents for Endo F3 digestions should be avoided.

The enzyme has a molecular mass of ~30 kDa.

The enzyme is provided as a solution in 20 mM Tris-HCl, pH 7.5.

Deglycosylation Procedure

1. Add up to 200 μg of glycoprotein to a microcentrifuge tube. Adjust the final volume to 37.5 μl with deionized water.
2. Add 10 μl of 5 \times Reaction Buffer (Catalog Number R9150)
3. Add 2.0 μl of Endo F3 solution to the reaction. Incubate 1 hour at 37°C.

Note: When cleaving non-core fucosylated biantennary and triantennary N-linked oligosaccharides, increase incubation time to 4 days. Many glycoproteins are contaminated with proteases. It is advised to add suitable protease inhibitors if long duration incubations are anticipated.

4. Monitor cleavage by SDS-PAGE.

Figure 1.
Specificities of Endoglycosidases F1, F2, and F3

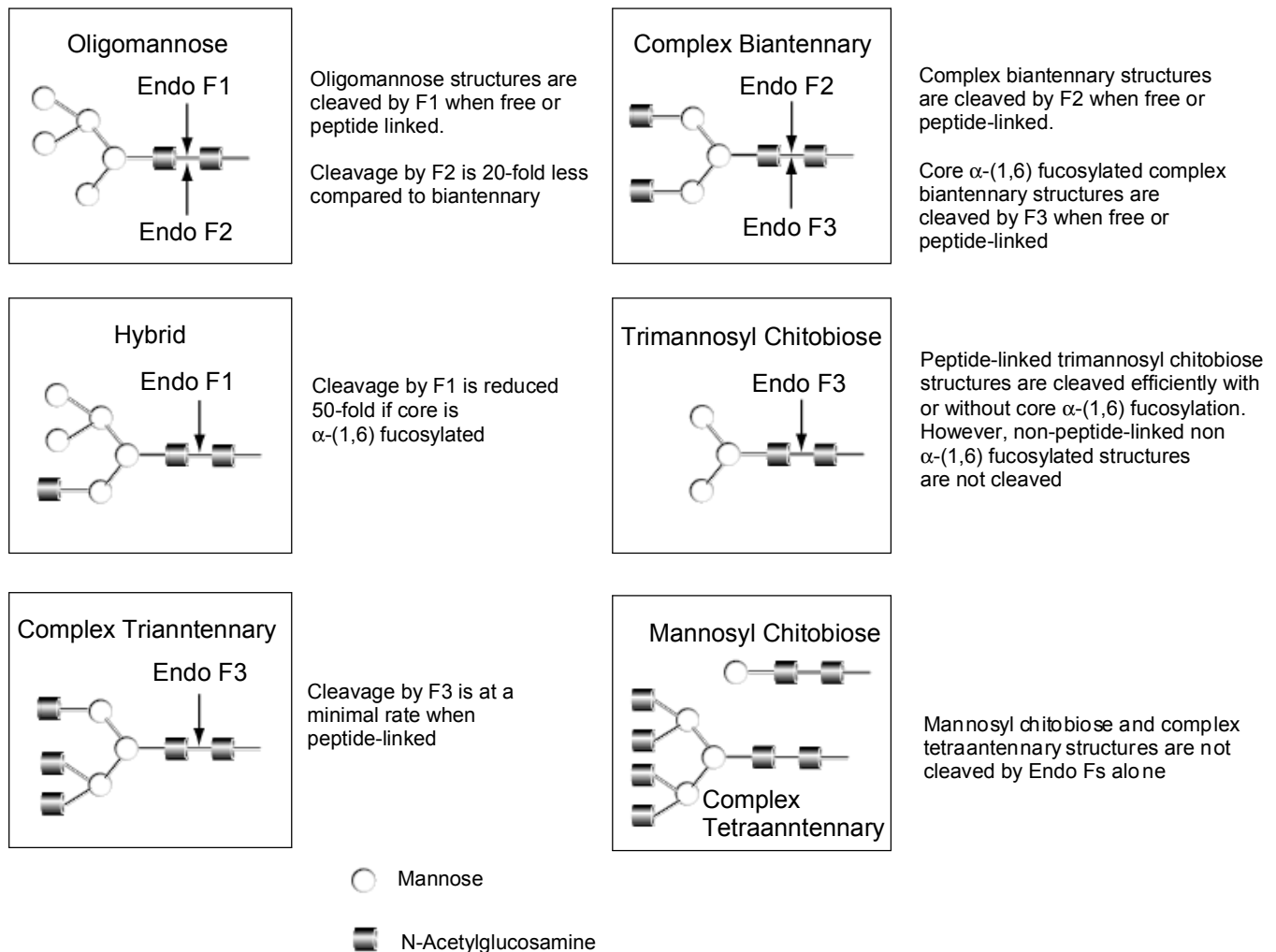


Figure 2.
Specificities of Endoglycosidases F1, F2, and F3 in conjunction with the EDEGLY Kit

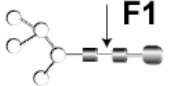
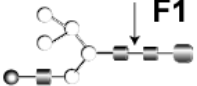
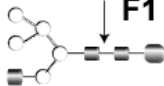
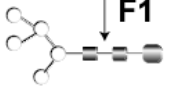
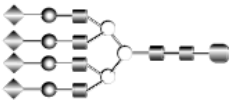
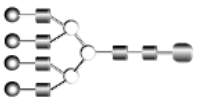
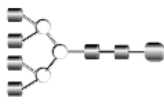
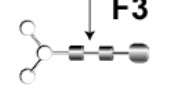
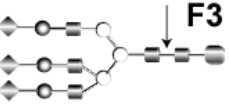
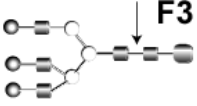
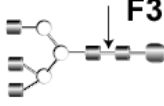

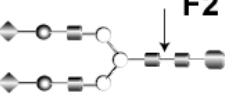
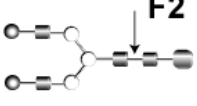
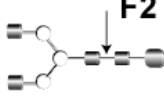






		With neuraminidase N8271	With β -galactosidase G0413	With β -N-acetylglucosaminidase A6805		
Oligomannose						
Hybrid						
Tetraantennary						
Triantennary						
Biantennary						
		 N-Acetylneuraminic Acid	 Galactose	 Mannose	 N-Acetylglucosamine	 Asn

Figure 3.
Specificity of Engoglycosidase F1

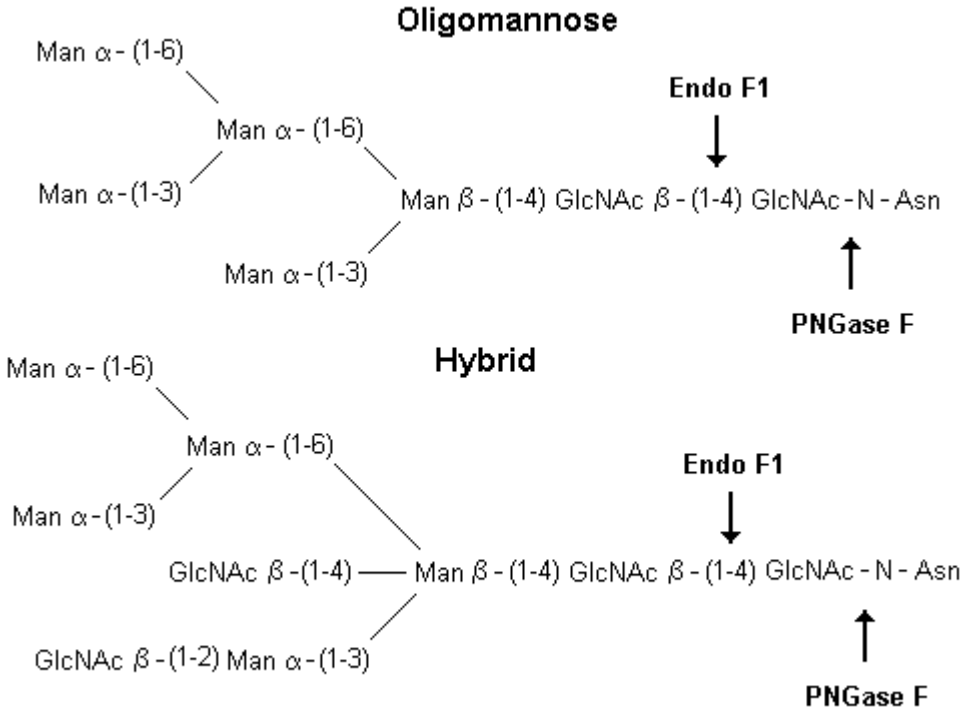
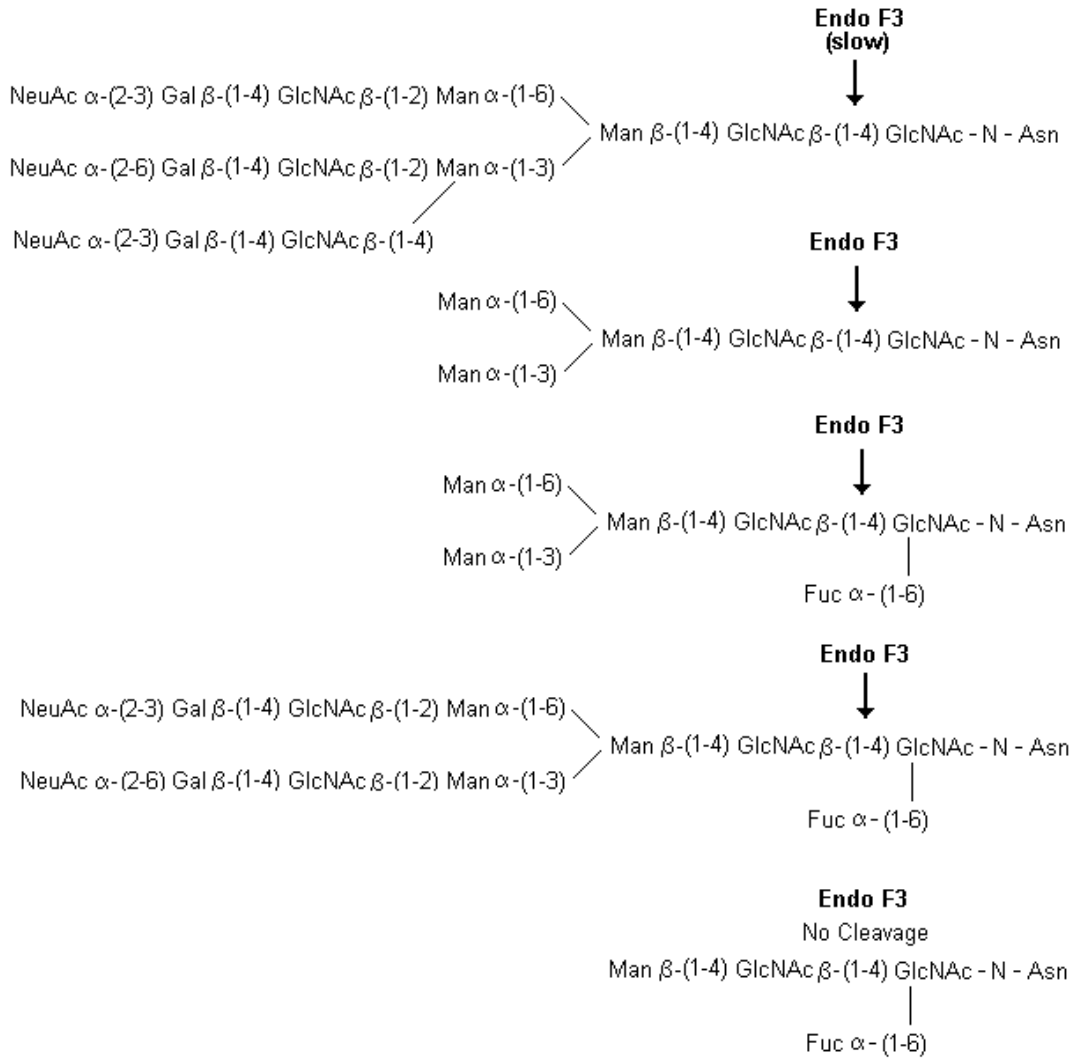


Figure 4.
Specificity of Endoglycosidase F3



References

1. Maley, P. *et al.*, Characterization of glycoproteins and their associated oligosaccharides through the use of endoglycosidases. *Anal. Biochem.*, **180**, 195-204 (1989).
2. Plummer, T.H. Jr. *et al.*, Porcine fibrinogen glycopeptides: substrates for detecting endo- β -N-cetylglucosaminidases F2 and F3. *Anal. Biochem.*, **235**, 98-101 (1996).
3. Reddy, A. *et al.*, High-level expression of the Endo-beta-N-acetylglucosaminidase F2 gene in *E. coli*: one step purification to homogeneity. *Glycobiology*, **8**, 633-636 (1998).
4. Tarentino, A.L. *et al.*, Deglycosylation of Asparagine-Linked Glycans by Peptide: N-Glycosidase F. *Biochemistry*, **24**, 4665-4671 (1985).
5. Tarentino, A.L. *et al.*, Multiple endoglycosidase (Endo) F activities expressed by *Flavobacterium meningosepticum*. Endo F1: molecular cloning, primary sequence, and structural relationship to Endo H. *J. Biol. Chem.*, **267**, 3868-3872 (1992).
6. Tarentino, A.L. *et al.*, Multiple endoglycosidase F activities expressed by *Flavobacterium meningosepticum* endoglycosidases F2 and F3: Molecular cloning, primary sequence, and enzyme expression. *J. Biol. Chem.*, **268**, 9702-9708 (1993).
7. Tarentino, A.L., and Plummer, T.H. Jr., Substrate specificity of *Flavobacterium meningosepticum*: Endo F2 and endo F3: purity is the name of the game. *Glycobiology*, **4**, 771-773 (1994).
8. Tarentino, A.L. and Plummer, T.H. Jr., Enzymatic deglycosylation of asparagine-linked glycans: purification, properties and specificity of oligosaccharide-cleaving enzymes from *Flavobacterium meningosepticum*. *Meth. Enzymol.*, **230**, 44-57 (1994).
9. Tarentino A.L. *et al.*, Overexpression and purification of non-glycosylated recombinant endo- β -N-acetylglucosaminidase F3. *Glycobiology*, **5**, 599-601 (1995).
10. Trimble, R.B., and Tarentino, A.L., Identification of Distinct Endoglycosidase (Endo) Activities in *Flavobacterium meningosepticum*: Endo F1, Endo F2 and Endo F3. *J. Biol. Chem.*, **266**, 1646-1651 (1991).

RBG,MAM 01/11-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.