

# PRODUCT INFORMATION



## 1F-Fructofuranosylnystose

Item No. 34292

**CAS Registry No.:** 59432-60-9  
**Formal Name:** O-β-D-fructofuranosyl-(2→1)-  
O-β-D-fructofuranosyl-(2→1)-  
O-β-D-fructofuranosyl-(2→1)-  
β-D-fructofuranosyl, α-D-  
glucopyranoside

**Synonyms:** GF<sub>4</sub>, 1,1,1-Kestopentaose

**MF:** C<sub>30</sub>H<sub>52</sub>O<sub>26</sub>

**FW:** 828.7

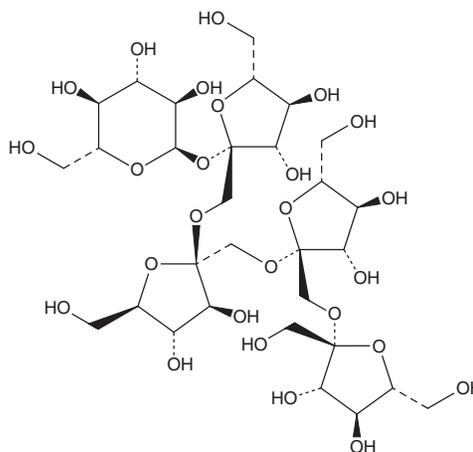
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Plant/*Morinda officinalis*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1F-Fructofuranosylnystose is supplied as a solid. A stock solution may be made by dissolving the 1F-fructofuranosylnystose in the solvent of choice, which should be purged with an inert gas. 1F-Fructofuranosylnystose is slightly soluble in DMSO and dimethyl formamide.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1F-fructofuranosylnystose can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1F-fructofuranosylnystose in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

1F-Fructofuranosylnystose is a fructooligosaccharide that has been found in onions.<sup>1</sup>

### Reference

1. Jaime, L., Martínez, F., Martín-Cabrejas, M.A., *et al.* Study of total fructan and fructooligosaccharide content in different onion tissues. *J. Sci. Food Agr.* **81(2)**, 177-182 (2000).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/30/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM