

# PRODUCT INFORMATION



## Xylotriose

Item No. 34033

CAS Registry No.: 47592-59-6

Formal Name: O-β-D-xylopyranosyl-(1→4)-O-β-D-xylopyranosyl-(1→4)-D-xylose

MF: C<sub>15</sub>H<sub>26</sub>O<sub>13</sub>

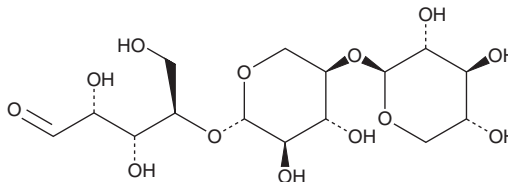
FW: 414.4

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

### Laboratory Procedures

Xylotriose is supplied as a solid. A stock solution may be made by dissolving the xylotriose in the solvent of choice, which should be purged with an inert gas. Xylotriose is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of xylotriose in these solvents is approximately 30 mg/ml.

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 xylotriose can be prepared by directly dissolving the solid in aqueous buffers. The solubility of xylotriose in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Xylotriose is a xylooligosaccharide that has been found in corn and various agricultural wastes.<sup>1,2</sup>

### References

1. Chapla, D., Pandit, P., and Shah, A. Production of xylooligosaccharides from corncob xylan by fungal xylanase and their utilization by probiotics. *Bioresour. Technol.* **115**, 215-221 (2012).
2. Akpinar, O., Erdogan, K., and Bostanci, S. Enzymatic production of Xylooligosaccharide from selected agricultural wastes. *Food Bioprod. Process.* **87(2)**, 145-151 (2009).

#### 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, 11/04/2022

#### 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