

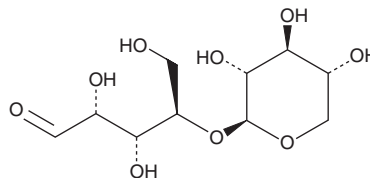
PRODUCT INFORMATION



1,4-D-Xylobiose

Item No. 34031

CAS Registry No.: 6860-47-5
Formal Name: 4-O-β-D-xylopyranosyl-D-xylose
Synonyms: 1,4-β-Xylobiose, β-(1,4)-Xylobiose
MF: C₁₀H₁₈O₉
FW: 282.2
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

1,4-D-Xylobiose is supplied as a solid. Aqueous solutions of 1,4-D-xylobiose can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1,4-D-xylobiose in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1,4-D-Xylobiose is a 1,4-linked disaccharide. It is a degradation product of hemicellulose, which is catabolized by xylanases.¹ It has been found in *T. reesei* endoxylanase-treated birch kraft pulp.²

References

1. Garg, S. Xylanase: Applications in biofuel production. *Curr. Metabolomics* **4**(1), 23-37 (2016).
2. Rydlund, A. and Dahlman, O. Oligosaccharides obtained by enzymatic hydrolysis of birch kraft pulp xylan: Analysis by capillary zone electrophoresis and mass spectrometry. *Carbohydr. Res.* **300**(2), 95-102 (1997).

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, 11/15/2022

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