

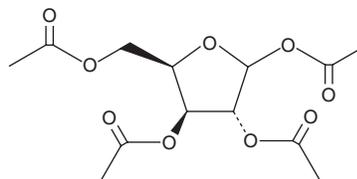
PRODUCT INFORMATION



1,2,3,5-Tetra-O-acetyl-D-xylofuranose

Item No. 17925

CAS Registry No.: 42927-46-8
Formal Name: 1,2,3,5-tetraacetate D-xylofuranose
MF: $C_{13}H_{18}O_9$
FW: 318.3
Purity: $\geq 95\%$
Supplied as: A neat oil
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

1,2,3,5-Tetra-O-acetyl-D-xylofuranose is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2,3,5-tetra-O-acetyl-D-xylofuranose in the solvent of choice, which should be purged with an inert gas. 1,2,3,5-Tetra-O-acetyl-D-xylofuranose is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1,2,3,5-tetra-O-acetyl-D-xylofuranose in ethanol and DMSO is approximately 15 mg/ml and approximately 30 mg/ml in DMF.

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 1,2,3,5-tetra-O-acetyl-D-xylofuranose can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 1,2,3,5-tetra-O-acetyl-D-xylofuranose in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1,2,3,5-Tetra-O-acetyl-D-xylofuranose is a starting material for the synthesis of nucleosides.¹

Reference

1. Singh, S.K., Sharma, R.K., Maity, J., *et al.* Greener biocatalytic approach to the synthesis of nucleosides and their precursors. *Nucleic Acids Symp. Ser. (Oxf.)* **52**, 273-274 (2008).

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, 11/23/2022

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