

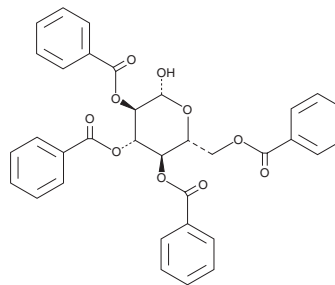
PRODUCT INFORMATION



2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose

Item No. 17163

CAS Registry No.: 64768-20-3
Formal Name: 2,3,4,6-tetrabenzoate β-D-glucopyranose
MF: C₃₄H₂₈O₁₀
FW: 596.6
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 274 nm
Supplied as: A crystalline 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

2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,3,4,6-tetra-O-benzoyl-β-D-glucopyranose in the solvent of choice, which should be purged with an inert gas. 2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2,3,4,6-tetra-O-benzoyl-β-D-glucopyranose in these solvents is approximately 30 mg/ml.

2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2,3,4,6-tetra-O-benzoyl-β-D-glucopyranose should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose has a solubility of approximately 0.33 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose is a D-glucopyranose derivative that has been used in glucosylation reactions.¹⁻⁴

References

1. Allevi, P., Anastasia, M., Ciuffreda, P., *et al.* The first direct method for C-glucopyranosyl derivatization of 2,3,4,6-tetra-O-benzyl-D-glucopyranose. *J. Chem. Soc. Chem. Commun.* **16**, 1245-1246 (1987).
2. Allevi, P., Anastasia, M., Ciuffreda, P., *et al.* C-Glucopyranosyl derivatives from readily available 2,3,4,6-tetra-O-benzyl-α-D-glucopyranosyl chloride. *J. Chem. Soc. Chem. Commun.* **2**, 101-102 (1987).
3. Koto, S., Sato, T., Morishima, N., *et al.* The glucosylation of several alcohols with tetra-O-benzyl-α-D-gluco-pyranose and a mixture of *p*-nitrobenzenesulfonyl chloride, silver triflorormethanesulfonate, and triethylamine. *Bull. Chem. Soc. Jpn.* **53**, 1761-1762 (1980).
4. Pavia, A.A., Rocheville, J.-M., and Ung, S.N. Nouvelle méthode de synthèse Stéréosélective de glycosides. Synthèse des α,α-tréhalose, analogues *galacto*, *manno* et autres α-D-glycosides. *Carb. Res.* **79(1)**, 79-89 (1980).

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, 11/10/2022

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