

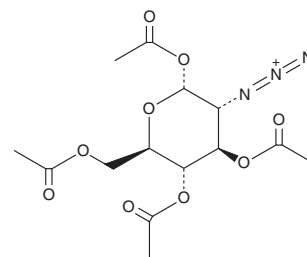
PRODUCT INFORMATION



1,3,4,6-Tetra-O-acetyl-2-azido-2-deoxy- α -D-Glucopyranose

Item No. 31153

CAS Registry No.: 56883-33-1
Formal Name: 2-azido-2-deoxy-1,3,4,6-tetraacetate- α -D-glucopyranose
MF: C₁₄H₁₉N₃O₉
FW: 373.3
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

1,3,4,6-Tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose is supplied as a solid. A stock solution may be made by dissolving the 1,3,4,6-tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose in the solvent of choice, which should be purged with an inert gas. 1,3,4,6-Tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 1,3,4,6-tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose 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 1,3,4,6-tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1,3,4,6-tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1,3,4,6-Tetra-O-acetyl-2-azido-2-deoxy- α -D-glucopyranose is a carbohydrate building block.¹ It has been used in the synthesis of amino-deoxy oligosaccharides.

Reference

1. Vasella, A., Witzig, C., Chiara, J.-L., *et al.* Convenient synthesis of 2-azido-2-deoxy-aldoses by diazo transfer. *Helv. Chim. Acta.* **74**(8), 2073-2077 (1991).

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