

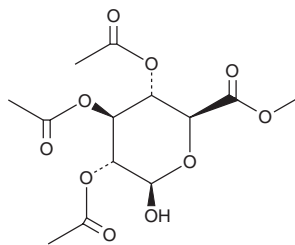
PRODUCT INFORMATION



2,3,4-Tri-O-acetyl-β-D-Glucuronide methyl ester

Item No. 30540

CAS Registry No.: 73464-50-3
Formal Name: 2,3,4-triacetate, β-D-glucopyranuronic acid, methyl ester
MF: C₁₃H₁₈O₁₀
FW: 334.3
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

2,3,4-Tri-O-acetyl-β-D-glucuronide methyl ester is supplied as a solid. A stock solution may be made by dissolving the 2,3,4-tri-O-acetyl-β-D-glucuronide methyl ester in the solvent of choice, which should be purged with an inert gas. 2,3,4-Tri-O-acetyl-β-D-glucuronide methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2,3,4-tri-O-acetyl-β-D-glucuronide methyl ester in ethanol and DMF is approximately 15 mg/ml and approximately 10 mg/ml in DMSO.

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 2,3,4-tri-O-acetyl-β-D-glucuronide methyl ester can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 2,3,4-tri-O-acetyl-β-D-glucuronide methyl ester in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2,3,4-Tri-O-acetyl-β-D-glucuronide methyl ester is a synthetic intermediate in the synthesis of prodrug forms of the antitumor antibiotic daunomycin (daunorubicin; Item No. 14159).¹

Reference

1. Leenders, R.G.G. and Scheeren, H.W. Synthesis and evaluation of novel daunomycin-phosphate-sulfate-β-glucuronide and -β-glucoside prodrugs for application in ADEPT. *Bioorg. Med. Chem. Lett.* **5**(24), 2975-2980 (1995).

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, 12/08/2022

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