

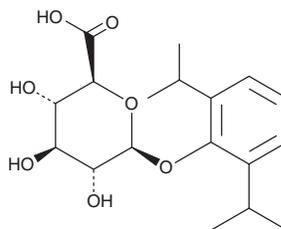
PRODUCT INFORMATION



Propofol β -D-Glucuronide

Item No. 19517

CAS Registry No.: 114991-26-3
Formal Name: 2,6-bis(1-methylethyl)phenyl β -D-glucopyranosiduronic acid
MF: C₁₈H₂₆O₇
FW: 354.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 211 nm
Supplied as: A crystalline 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

Propofol β -D-glucuronide is supplied as a crystalline solid. A stock solution may be made by dissolving the propofol β -D-glucuronide in the solvent of choice, which should be purged with an inert gas. Propofol β -D-glucuronide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of propofol β -D-glucuronide 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 propofol β -D-glucuronide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of propofol β -D-glucuronide in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Propofol β -D-glucuronide is a metabolite of the anesthetic and sedative propofol.¹ It is formed from propofol from the UDP-glucuronosyltransferase (UGT) isoform UGT1A9.

Reference

1. Chen, A., Zhou, X., Cheng, Y., *et al.* Design and optimization of the cocktail assay for rapid assessment of the activity of UGT enzymes in human and rat liver microsomes. *Toxicol. Lett.* **295**, 379-389 (2018).

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, 04/17/2023

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