

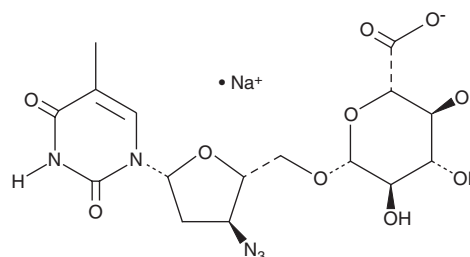
PRODUCT INFORMATION



3'-Azido-3'-deoxythymidine β -D-glucuronide (sodium salt)

Item No. 18679

CAS Registry No.: 133525-01-6
Formal Name: 3'-azido-3'-deoxy-5'-O- β -D-glucopyranuronosyl-thymidine, monosodium salt
Synonyms: AZT glucuronide, Zidovudine glucuronide
MF: C₁₆H₂₀N₅O₁₀ • Na
FW: 465.4
Purity: \geq 95%
UV/Vis.: λ_{max} : 267 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

3'-Azido-3'-deoxythymidine β -D-glucuronide (AZT glucuronide) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the AZT glucuronide (sodium salt) in the solvent of choice, which should be purged with an inert gas. AZT glucuronide (sodium salt) is soluble in the organic solvent DMSO at a concentration of approximately 0.5 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 AZT glucuronide (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of AZT glucuronide (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AZT glucuronide is a metabolite of the nucleoside reverse transcriptase inhibitor zidovudine (Item No. 15492).¹ It is formed from zidovudine by the UDP-glucuronosyltransferase (UGT) isoform UGT2B7.

Reference

1. Walsky, R.L., Bauman, J.N., Bourcier, K., *et al.* Optimized assays for human UDP-glucuronosyltransferase (UGT) activities: Altered alamethicin concentration and utility to screen for UGT inhibitors. *Drug Metab. Dispos.* **40**(5), 1051-1065 (2012).

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, 06/23/2023

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