

PRODUCT INFORMATION



UDP-Glucuronic Acid (sodium salt hydrate)

Item No. 20674

Formal Name: α -D-glucopyranuronic acid, 1 \rightarrow P'-ester with uridine 5'-(trihydrogen diphosphate), trisodium salt, hydrate

Synonyms: UDPGA, UDP-GlcA, Uridine-5'-diphosphoglucuronic Acid

MF: C₁₅H₁₉N₂O₁₈P₂ • 3Na [XH₂O]

FW: 646.2

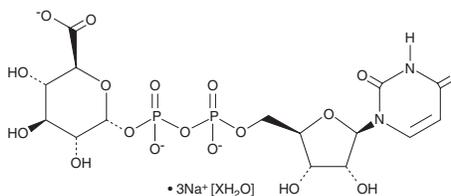
Purity: \geq 95%

UV/Vis.: λ_{max} : 262 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

UDP-Glucuronic acid (sodium salt hydrate) is supplied as a crystalline solid. Aqueous solutions of UDP-glucuronic acid (sodium salt hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of UDP-glucuronic acid (sodium salt hydrate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

UDP-Glucuronic acid is a nucleoside diphosphate sugar that is a source of glucuronic acid for polysaccharide biosynthesis. This compound can be used for the formation of many glucuronides with various aglycones.¹

Reference

1. Green, M.D. and Tephly, T.R. Glucuronidation of amine substrates by purified and expressed UDP-glucuronosyltransferase proteins. *Drug Metab. Dispos.* **26(9)**, 860-867 (1998).

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, 11/18/2022

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