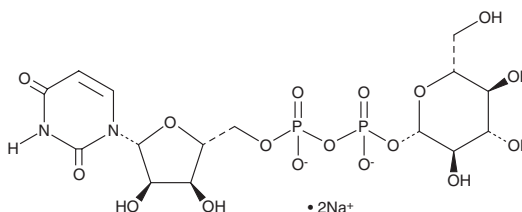


PRODUCT INFORMATION

UDP-β-D-Glucose (sodium salt)

Item No. 21620

CAS Registry No.: 7333-33-7
Formal Name: uridine 5'-(trihydrogen pyrophosphate), mono-β-D-glucopyranosyl ester, disodium salt
Synonyms: NSC 20269, UDPG, UDP-Glucose, Uridine-5'-diphosphoglucose
MF: C₁₅H₂₂N₂O₁₇P₂ • 2Na
FW: 610.3
Purity: ≥95%
UV/Vis.: λ_{max}: 261 nm
Supplied as: A crystalline 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

UDP-β-D-Glucose (sodium salt) is supplied as a crystalline solid. Aqueous solutions of UDP-β-D-glucose (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of UDP-β-D-glucose (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

UDP-β-D-glucose is a nucleotide sugar, purinergic P2Y₁₄ receptor agonist, and β-anomer of UDP-α-D-glucose (Item No. 15602).¹ It induces inositol phosphate production in COS-7 cells expressing the human P2Y₁₄ receptor (EC₅₀ = 0.588 μM).

Reference

1. Ko, H., Das, A., Carter, R.L., *et al.* Molecular recognition in the P2Y₁₄ receptor: Probing the structurally permissive terminal sugar moiety of uridine-5'-diphosphoglucose. *Bioorg. Med. Chem.* **17(14)**, 5298-5311 (2009).

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, 04/03/2025

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