

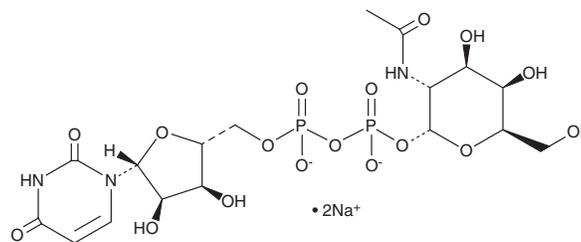
PRODUCT INFORMATION



UDP-GalNAc (sodium salt)

Item No. 36506

CAS Registry No.: 108320-87-2
Formal Name: uridine 5'-(trihydrogen diphosphate), P'-[2-(acetylamino)-2-deoxy- α -D-galactopyranosyl] ester, disodium salt
Synonyms: UDP-N-acetyl-D-Galactosamine, Uridine 5'-diphosphate GalNAc
MF: $C_{17}H_{25}N_3O_{17}P_2 \cdot 2Na$
FW: 651.3
Purity: $\geq 90\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

UDP-GalNAc (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the UDP-GalNAc (sodium salt) in water. The solubility of UDP-GalNAc (sodium salt) in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

UDP-GalNAc is a nucleotide sugar that is used for glycosylation by glycosyltransferases, which transfer the N-acetyl-D-galactosamine (Item No. 31728) moiety to various substrates.^{1,2}

References

1. Yazer, M.H., Olsson, M.L., and Palcic, M.M. The *cis*-AB blood group phenotype: Fundamental lessons in glycobiology. *Transfus. Med. Rev.* **20(3)**, 207-217 (2006).
2. Wang, Y., Abernethy, J.L., Eckhardt, A.E., *et al.* Purification and characterization of a UDP-GalNAc:polypeptide N-acetylgalactosaminyltransferase specific for glycosylation of threonine residues. *J. Biol. Chem.* **267(18)**, 12709-12716 (1992).

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, 10/21/2025

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