

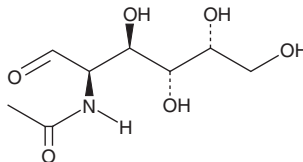
PRODUCT INFORMATION



N-acetyl-D-Galactosamine

Item No. 31728

CAS Registry No.: 1811-31-0
Formal Name: 2-(acetylamino)-2-deoxy-D-galactose
Synonyms: GalNAc, N-Acetylgalactosamine
MF: C₈H₁₅NO₆
FW: 221.2
Purity: ≥95%
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

N-acetyl-D-Galactosamine is supplied as a crystalline solid. Aqueous solutions of N-acetyl-D-galactosamine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N-acetyl-D-galactosamine in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-acetyl-D-Galactosamine (GalNAc) is an amino sugar derivative of galactose and a component of O-GalNAc glycans.^{1,2} It is transferred from UDP-GalNAc to the hydroxy group of protein serine or threonine residues by polypeptide GalNAc transferase, forming the Tn antigen, during the first step of mucin-type O-glycosylation.^{2,3} GalNAc has been used as a targeting moiety for liver-targeted delivery of oligonucleotides.¹

References

1. Huang, Y. Preclinical and clinical advances of GalNAc-decorated nucleic acid therapeutics. *Mol. Ther. Nucleic Acids* **6**, 116-132 (2017).
2. Brockhausen, I., Schachter, H., and Stanley, P. O-GalNAc glycans. *Essentials of Glycobiology - NCBI Bookshelf*. Varki, A., Cummings, R.D., Esko, J.D., et al. editors, 2nd edition, Harbor Laboratory Press (2009).
3. Bennett, E.P., Mandel, U., Clausen, H., et al. Control of mucin-type O-glycosylation: A classification of the polypeptide GalNAc-transferase gene family. *Glycobiology* **22(6)**, 736-756 (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, 11/02/2022

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