

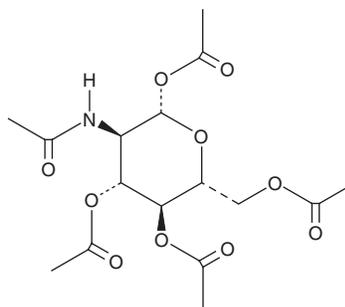
PRODUCT INFORMATION



β-D-Glucosamine Pentaacetate

Item No. 17867

CAS Registry No.: 7772-79-4
Formal Name: 2-(acetylamino)-2-deoxy-β-D-glucopyranose, 1,3,4,6-tetraacetate
Synonyms: N-Acetyl-β-D-Glucosamine tetraacetate, NSC 224432
MF: C₁₆H₂₃NO₁₀
FW: 389.4
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

β-D-Glucosamine pentaacetate is supplied as a crystalline solid. A stock solution may be made by dissolving the β-D-glucosamine pentaacetate in the solvent of choice. β-D-Glucosamine pentaacetate is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of β-D-glucosamine pentaacetate in these solvents is approximately 20 and 30 mg/ml, respectively.

β-D-Glucosamine pentaacetate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, β-D-glucosamine pentaacetate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. β-D-Glucosamine pentaacetate has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

β-D-Glucosamine pentaacetate is an N-acetylglucosamine derivative that has been shown to promote hyaluronic acid production.^{1,2}

References

1. Bastida, A., Sabuquillo, P., Armisen, P., *et al.* A single step purification, immobilization, and hyperactivation of lipases *via* interfacial adsorption on strongly hydrophobic supports. *Biotechnol. Bioeng.* **58(5)**, 486-493 (1998).
2. Schulz, T., Schumacher, U., and Prehm, P. Hyaluronan export by the ABC transporter MRP5 and its modulation by intracellular cGMP. *J. Biol. Chem.* **282(29)**, 20999-21004 (2007).

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, 12/14/2022

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