

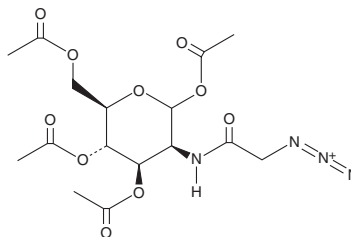
PRODUCT INFORMATION



Ac4ManNAz

Item No. 36283

CAS Registry No.: 361154-30-5
Formal Name: 2-[(2-azidoacetyl)amino]-2-deoxy-D-mannopyranose, 1,3,4,6-tetraacetate
Synonyms: N-Azidoacetyl-D-mannosamine, N-Azidoacetylmannosamine-tetraacetylated
MF: C₁₆H₂₂N₄O₁₀
FW: 430.4
Purity: ≥98%
Supplied as: A 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

Ac4ManNAz is supplied as a solid. A stock solution may be made by dissolving the Ac4ManNAz in the solvent of choice, which should be purged with an inert gas. Ac4ManNAz is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Ac4ManNAz in these solvents is approximately 2 mg/ml. Ac4ManNAz is also slightly soluble in ethanol.

Description

Ac4ManNAz is an azide-functionalized monosaccharide.^{1,2} It has been used for metabolic labeling to study the cellular trafficking of labeled proteins.

References

1. Han, S.-S., Shim, H.-E., Park, S.-J., *et al.* Safety and optimization of metabolic labeling of endothelial progenitor cells for tracking. *Sci. Rep.* **8**(1), 13212 (2018).
2. Rochefort, M., Girgis, M.D., Ankeny, J.S., *et al.* Metabolic exploitation of the sialic acid biosynthetic pathway to generate site-specifically labeled antibodies. *Glycobiology* **24**(1), 62-69 (2014).

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, 12/19/2022

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