

PRODUCT INFORMATION



Chitobiose Octaacetate

Item No. 16379

CAS Registry No.: 41670-99-9
Formal Name: 2-(acetylamino)-2-deoxy-4-O-[3,4,6-tri-O-acetyl-2-(acetylamino)-2-deoxy-β-D-glucopyranosyl]-D-glucopyranose-1,3,6-triacetate

Synonym: Peracetylchitobiose

MF: C₂₈H₄₀N₂O₁₇

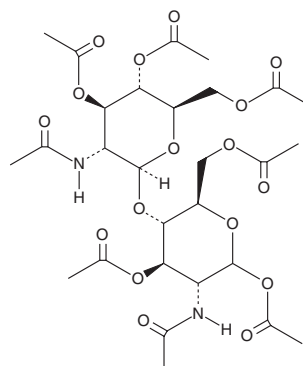
FW: 676.6

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

Chitobiose octaacetate is supplied as a crystalline solid. A stock solution may be made by dissolving the chitobiose octaacetate in the solvent of choice, which should be purged with an inert gas. Chitobiose octaacetate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of chitobiose octaacetate in ethanol is approximately 0.5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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

Description

Chitobiose is a dimer of β-1,4 linked glucosamine units derived from chitin, which, in turn, is a long chain polymer of N-acetylglucosamine that is a primary component of fungal cell walls and arthropod exoskeletons. Chitobiose octaacetate is a form of chitobiose containing eight acetate groups. It is formed by the acetolysis of chitin. Chitobiose octaacetate can be used for the synthesis of chitobiose oxazoline and other chitin precursors.

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