

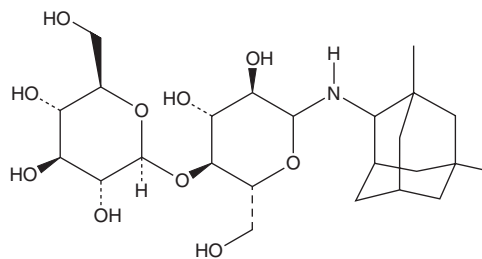
PRODUCT INFORMATION



Memantine Lactose Adduct

Item No. 20662

CAS Registry No.: 1159637-28-1
Formal Name: N-(3,5-dimethyltricyclo[3.3.1.1^{3,7}]dec-1-yl)-4-O-β-D-glucopyranosyl-D-glucopyranosylamine
MF: C₂₄H₄₁NO₁₀
FW: 503.6
Purity: ≥95%
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

Memantine lactose adduct is supplied as a solid. A stock solution may be made by dissolving the memantine lactose adduct in the solvent of choice, which should be purged with an inert gas. Memantine lactose adduct is soluble in the organic solvent methanol. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Memantine lactose adduct is a lactose adduct of memantine (Item No. 14184) and an intermediate in the synthesis of adamantane amine antibiotics.¹

Reference

1. Assenza, S.P., Shen, N., Iacono, S., et al. Reducing carbohydrate derivatives of adamantane amines, and synthesis and methods of use thereof. **US 20060211650 A1**, (2006).

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, 12/16/2022

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