

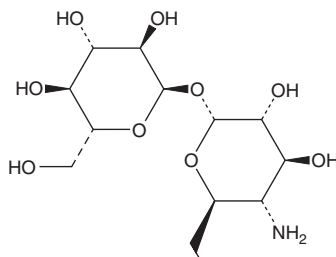
PRODUCT INFORMATION



4-Trehalosamine

Item No. 40229

CAS Registry No.: 51855-99-3
Formal Name: α -D-glucopyranosyl 4-amino-4-deoxy- α -D-glucopyranoside
MF: $C_{12}H_{23}NO_{10}$
FW: 341.3
Purity: $\geq 90\%$
Supplied as: A powder
Storage: $-20^{\circ}C$
Stability: ≥ 1 year
Item Origin: Bacterium/*Streptomyces* sp.



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

4-Trehalosamine is supplied as a powder. A stock solution may be made by dissolving the 4-trehalosamine in the solvent of choice. 4-Trehalosamine is soluble in organic solvents such as DMSO and methanol, which should be purged with an inert gas. 4-Trehalosamine is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

4-Trehalosamine is a microbial metabolite that has been found in *Streptomyces* and has antibiotic activity.¹ It is active against *E. coli*, *K. pneumoniae*, and *B. subtilis* in cup assays when used at a concentration of 0.5, 1, and 2 mg/ml.

Reference

1. Naganawa, H., Usui, N., Takita, T., *et al.* 4-Amino-4-deoxy- α , α -trehalose, a new metabolite of a *Streptomyces*. *J. Antibiot. (Tokyo)* **27**(2), 145-146 (1974).

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, 11/07/2023

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