

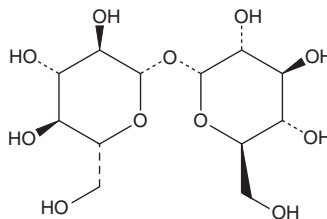
PRODUCT INFORMATION



α,β -Trehalose

Item No. 44333

CAS Registry No.: 585-91-1
Formal Name: β -D-glucopyranosyl α -D-glucopyranoside
MF: $C_{12}H_{22}O_{11}$
FW: 342.3
Purity: $\geq 95\%$
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

α,β -Trehalose is supplied as a solid. Aqueous solutions of α,β -trehalose can be prepared by directly dissolving the solid in aqueous buffers. α,β -Trehalose is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

α,β -Trehalose is a disaccharide that has been found in *C. cucullus*.¹

Reference

1. Sogame, Y., Ogata, M., Hakozaiki, S., *et al.* α,β -trehalose, an intracellular substance in resting cyst of colpodid ciliates as a key to environmental tolerances. *Biochem. Biophys. Res. Commun.* **716**, 149971 (2024).

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, 11/07/2025

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