

PRODUCT INFORMATION



Trehalose C14

Item No. 44264

CAS Registry No.: 64622-92-0

Formal Name: α -D-glucopyranoside, α -D-glucopyranosyl, 6-tetradecanoate

Synonym: Trehalose 6-myristate

MF: $C_{26}H_{48}O_{12}$

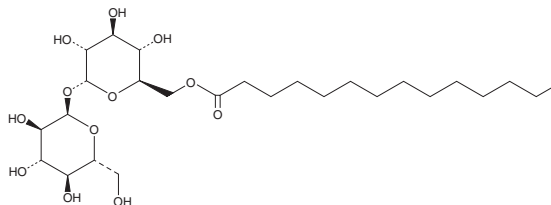
FW: 552.7

Purity: $\geq 98\%$

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

Trehalose C14 is supplied as a solid. A stock solution may be made by dissolving the trehalose C14 in the solvent of choice, which should be purged with an inert gas. Trehalose C14 is soluble (≥ 10 mg/ml) in DMSO. Trehalose C14 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

Trehalose C14 is a disaccharide-based surfactant that has been used in the solubilization of cell membranes for lipidomics analysis and as an emulsifier.^{1,2} It has a critical micelle concentration (CMC) of 0.012 mM.

References

1. Sakata, J., Furusho, A., Sugiyama, E., *et al.* Development of a highly efficient solubilization method for mass spectrometric analysis of phospholipids in living single cells. *Anal. Sci.* **40**(5), 917-924 (2024).
2. Nakayama, S., Kimura, Y., Miki, S., *et al.* Influence of sugar surfactant structure on the encapsulation of oil droplets in an amorphous sugar matrix during freeze-drying. *Food Res. Int.* **70**, 143-149 (2015).

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, 09/18/2025

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