

# PRODUCT INFORMATION



## Lentztrehalose C

Item No. 37734

**CAS Registry No.:** 1808096-68-5  
**Formal Name:**  $\alpha$ -D-glucopyranosyl 3,4-O-[(1R)-1-(1-hydroxy-1-methylethyl)-1,2-ethanediyl]- $\alpha$ -D-glucopyranoside

**MF:** C<sub>17</sub>H<sub>30</sub>O<sub>12</sub>

**FW:** 426.4

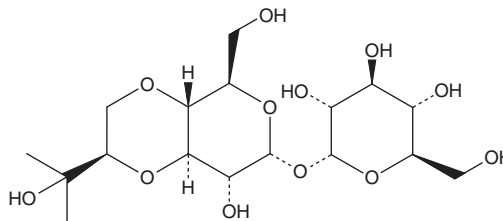
**Purity:**  $\geq 75\%$

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq 4$  years

**Item Origin:** Bacterium/*Lentzea* sp. ML457-mF



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

### Laboratory Procedures

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

### Description

Lentztrehalose C is a disaccharide that has been found in *Lentzea* and has autophagy-inducing activity.<sup>1</sup> It induces autophagy in MeWo melanoma and OVK18 ovarian cancer cells when used at a concentration of 100 mM.

### Reference

1. Wada, S., Kubota, Y., Sawa, R., *et al.* Novel autophagy inducers lentztrehaloses A, B and C. *J. Antibiot. (Tokyo)* **68(8)**, 521-529 (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, 06/13/2023

#### 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