

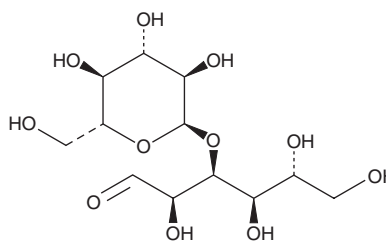
PRODUCT INFORMATION



Nigerose

Item No. 37419

CAS Registry No.: 497-48-3
Formal Name: 3-O- α -D-glucopyranosyl-D-glucose
Synonym: Sakebiose
MF: $C_{12}H_{22}O_{11}$
FW: 342.3
Purity: $\geq 90\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

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

Description

Nigerose is a disaccharide that has been found in *Acremonium*.¹ It prevents increases in the levels of malondialdehyde (MDA) and reactive oxygen species (ROS) and inhibits lactate dehydrogenase (LDH) release in HL-60 cells post-cryopreservation when used at a concentration of 150 μM in combination with the glycoside salidroside (Item No. 13628).²

References

1. Konishi, Y. and Shindo, K. Production of nigerose, nigerosyl glucose, and nigerosyl maltose by *Acremonium* sp. S4G13. *Biosci. Biotechnol. Biochem.* **61**(3), 439-442 (1997).
2. Straka, M.-S., Al-Otaibi, N.A., Whitfield, P.D., et al. Disaccharides-based cryo-formulant effect on modulating phospho/mitochondrial lipids and biological profiles of human leukaemia cells. *Cell. Physiol. Biochem.* **55**(2), 206-221 (2021).

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, 06/01/2023

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