

PRODUCT INFORMATION



Siamenoside I

Item No. 34656

CAS Registry No.: 126105-12-2
Formal Name: (3 β ,9 β ,10 α ,11 α ,24R)-3-(β -D-glucopyranosyloxy)-11,25-dihydroxy-9-methyl-19-norlanost-5-en-24-yl O- β -D-glucopyranosyl-(1 \rightarrow 2)-O-[β -D-glucopyranosyl-(1 \rightarrow 6)]- β -D-glucopyranoside

MF: C₅₄H₉₂O₂₄

FW: 1,125.3

Purity: \geq 90%

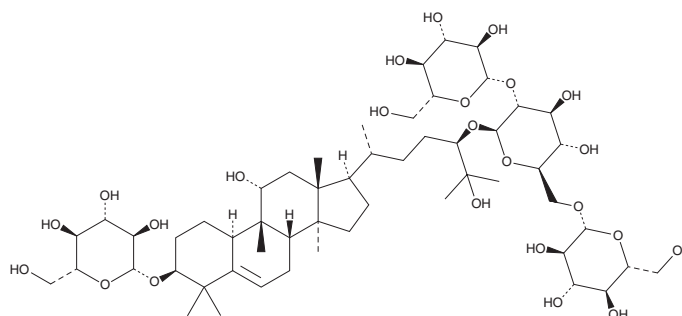
Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Momordica grosvenori*

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



Laboratory Procedures

Siamenoside I is supplied as a solid. A stock solution may be made by dissolving the siamenoside I in the solvent of choice, which should be purged with an inert gas. Siamenoside I is soluble in the organic solvent ethanol.

Description

Siamenoside I is a cucurbitane glycoside that has been found in *S. grosvenori* (monk fruit).¹ Formulations containing siamenoside I have been used as natural sweeteners.

Reference

1. Younes, M., Aquilina, G., Engel, K.-H., *et al.* Safety of use of monk fruit extract as a food additive in different food categories. *EFSA J.* **17(12)**, e05921 (2019).

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, 11/14/2022

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