

PRODUCT INFORMATION



11-oxo MGV

Item No. 22416

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

Synonym: 11-oxo Mogroside V

MF: C₆₀H₁₀₀O₂₉

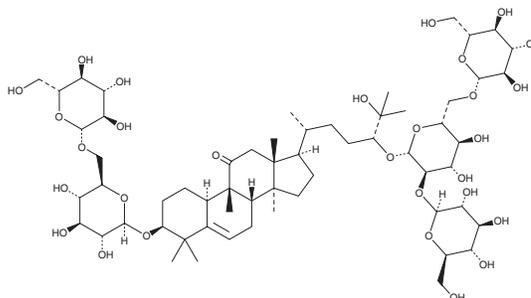
FW: 1,285.4

Purity: \geq 95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

11-oxo MGV is supplied as a crystalline solid. A stock solution may be made by dissolving the 11-oxo MGV in the solvent of choice, which should be purged with an inert gas. 11-oxo MGV is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 11-oxo MGV in these solvents is approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 11-oxo MGV can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 11-oxo MGV in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

11-oxo MGV is a sweet tasting cucurbitane glycoside derivative of mogroside V (Item No. 19853). It has antioxidant effects against both O₂⁻ and OH⁻ (EC₅₀s = 4.79 and 16.52 μ g/ml, respectively) and prevents OH-induced DNA damage (EC₅₀ = 3.09 μ g/ml).¹ 11-oxo MGV also inhibits carcinogenesis in murine models.²

References

1. Chen, W.J., Wang, J., and Qi, X.Y. The antioxidant activities of natural sweeteners, mogrosides, from fruits of *Siraitia grosvenori*. *Int. J. Food Sci. Nutr.* **58(7)**, 548-556 (2007).
2. Takasaki, M., Konoshima, T., Murata, Y., et al. Anticarcinogenic activity of natural sweeteners, cucurbitane glycosides, from *Momordica grosvenori*. *Cancer Lett.* **198(1)**, 37-42 (2003).

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, 02/08/2024

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