

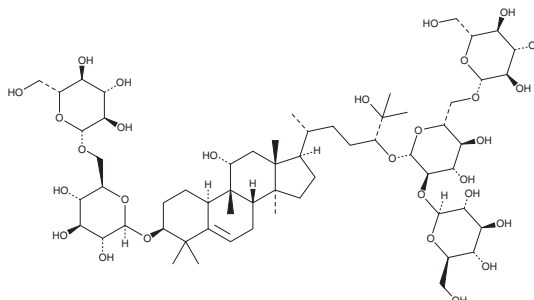
PRODUCT INFORMATION



Mogroside V Item No. 19853

CAS Registry No.: 88901-36-4
Formal Name: (3 β ,9 β ,10 α ,11 α ,24R)-3-[(6-O- β -D-glucopyranosyl- β -D-glucopyranosyl)oxy]-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

Synonym: MGV
MF: C₆₀H₁₀₂O₂₉
FW: 1,287.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 264 nm
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

Mogroside V is supplied as a crystalline solid. A stock solution may be made by dissolving the mogroside V in the solvent of choice, which should be purged with an inert gas. Mogroside V is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of mogroside V 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 mogroside V can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of mogroside V in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Mogroside V is a natural cucurbitane glycoside that can be isolated from *S. grosvenorii*.¹ It is a sweet tasting compound that is used as a low calorie sweetener.^{1,2} Mogroside V has antioxidant properties and possesses tumor growth inhibitory activity in pancreatic cancer models.^{2,3}

References

1. Zhang, M., Yang, H., Zhang, H., *et al.* Development of a process for separation of mogroside V from *Siraitia grosvenorii* by macroporous resins. *Molecules* **16**(9), 7288-7301 (2011).
2. Liu, C., Dai, L.-H., Dou, D.-Q., *et al.* A natural food sweetener with anti-pancreatic cancer properties. *Oncogenesis* **2015**(5) (2016).
3. Wang, L., Yang, Z., Lu, F., *et al.* Cucurbitane glycosides derived from mogroside II_E: Structure-taste relationships, antioxidant activity, and acute toxicity. *Molecules* **19**(8), 12676-12689 (2014).

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, 02/07/2024

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