

PRODUCT INFORMATION



Crocin II

Item No. 42485

CAS Registry No.: 55750-84-0

Formal Name: 1-[β-D-glucopyranosyl
(2E,4E,6E,8E,10E,12E,14E)-2,6,11,15-
tetramethyl-2,4,6,8,10,12,14-
hexadecaheptaenedioate]6-O-β-D-
glucopyranosyl-β-D-glucopyranose
Synonyms: *trans*-3-Gg-Crocin,
Crocin Gentiobiosylglucosyl ester,
Tricrocin

MF: C₃₈H₅₄O₁₉

FW: 814.8

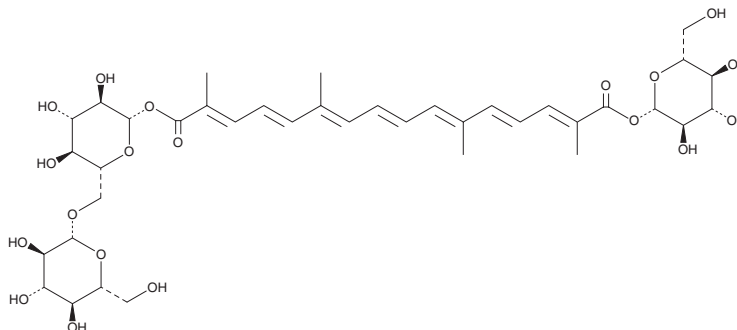
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: *Crocus sativus* L



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

Laboratory Procedures

Crocin II is supplied as a solid. A stock solution may be made by dissolving the crocin II in the solvent of choice, which should be purged with an inert gas. Crocin II is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Crocin II is a carotenoid glycoside that has been found in *G. jasminoides* and has antioxidant and proliferation-enhancing activities.^{1,2} It scavenges peroxy radicals in an oxygen radical scavenging capacity (ORAC) assay when used at concentrations of 10 and 20 μM.¹ Crocin II (5, 25, and 50 μg/ml) increases the proliferation rate of primary human mesenchymal stem cells (hMSCs).² It also reduces senescence in primary hMSCs when used at a concentration of 25 μg/ml.

References

1. Zhuang, G.-D., Gu, W.-T., Xu, S.-H., *et al.* Rapid screening of antioxidant from natural products by AAPH-Incubating HPLC-DAD-HR MS/MS method: A case study of *Gardenia jasminoides* fruit. *Food Chem.* **401**, 401:134091 (2023).
2. Yuan, T.-J., Xu, X.-H., Zhou, N., *et al.* Phytochemicals as new therapeutic candidates simultaneously stimulate proliferation and counteract senescence of stem cells. *Biomed Pharmacother.* **151**, 113170 (2022).

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, 12/17/2024

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