

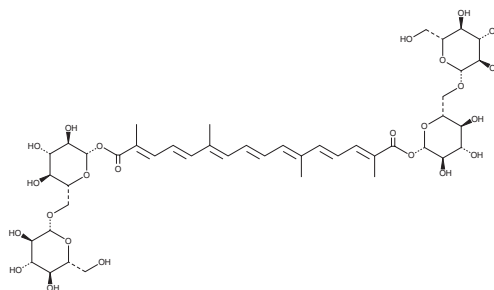
PRODUCT INFORMATION



Crocin/*Gardenia* Extract

Item No. 26337

CAS Registry No.: 42553-65-1
MF: C₄₄H₆₄O₂₄ (for crocin I)
FW: 977.0
UV/Vis.: λ_{max}: 324, 441 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Gardenia*



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

Laboratory Procedures

Crocin/*Gardenia* extract is supplied as a crystalline solid. A stock solution may be made by dissolving the crocin/*Gardenia* extract in the solvent of choice. Crocin/*Gardenia* extract is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of crocin/*Gardenia* extract in these solvents is approximately 10 and 5 mg/ml, respectively. Crocin/*Gardenia* extract is also partially soluble in ethanol.

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 crocin/*Gardenia* extract can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of crocin/*Gardenia* extract in PBS (pH 7.2) is approximately 3.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Crocin/*Gardenia* extract is a *Gardenia* extract that contains crocin I, II, and III, and crocetin, among other compounds.

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, 01/02/2025

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