

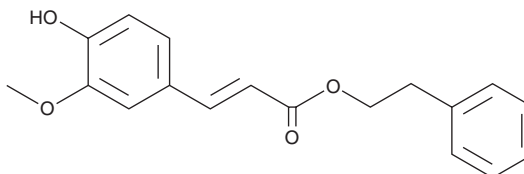
PRODUCT INFORMATION



Phenylethyl 3-methylcaffeate

Item No. 39152

CAS Registry No.: 71835-85-3
Formal Name: 3-(4-hydroxy-3-methoxyphenyl)-2-propenoic acid, 2-phenylethyl ester
Synonyms: NSC 666255, NSC 671193, Phenethyl Ferulate, 2-Phenylethyl Ferulate, β -Phenethyl Ferulate
MF: C₁₈H₁₈O₄
FW: 298.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Phenylethyl 3-methylcaffeate is supplied as a solid. A stock solution may be made by dissolving the phenylethyl 3-methylcaffeate in the solvent of choice, which should be purged with an inert gas. Phenylethyl 3-methylcaffeate is soluble in acetonitrile and methanol.

Description

Phenylethyl 3-methylcaffeate is a caffeic acid ester that has been found in propolis, a resin produced by honeybees, and has anticancer activity.^{1,2} Dietary administration of phenylethyl 3-methylcaffeate (750 ppm) reduces the incidence of both invasive and non-invasive colonic adenocarcinomas and the activity of colonic mucosal and tumor phosphatidylinositol-specific phospholipase C (PI-PLC) in a rat model of azoxymethane-induced carcinogenesis.¹ It also increases intratumoral apoptosis in a rat model of azoxymethane-induced carcinogenesis.²

References

1. Rao, C.V., Desai, D., Rivenson, A., *et al.* Chemoprevention of colon carcinogenesis by phenylethyl-3-methylcaffeate. *Cancer Res.* **55**(11), 2310-2315 (1995).
2. Samaha, H.S., Kelloff, G.J., Steele, V., *et al.* Modulation of apoptosis by sulindac, curcumin, phenylethyl-3-methylcaffeate, and 6-phenylhexyl isothiocyanate: apoptotic index as a biomarker in colon cancer chemoprevention and promotion. *Cancer Res.* **57**(7), 1301-1305 (1997).

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, 03/19/2024

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