

PRODUCT INFORMATION



Caffeic Acid phenylethyl ester

Item No. 70750

CAS Registry No.: 104594-70-9
Formal Name: (E)-3-(3,4-dihydroxyphenyl)-2-propenoic acid, 2-phenylethyl ester

Synonyms: CAPE, 2-Phenylethyl Caffeate, β -Phenylethyl Caffeate

MF: C₁₇H₁₆O₄

FW: 284.3

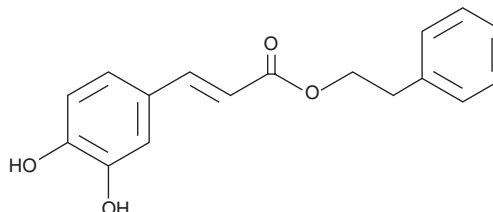
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 246, 300, 331 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

Caffeic acid phenylethyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the caffeic acid phenylethyl ester in the solvent of choice, which should be purged with an inert gas. Caffeic Acid phenylethyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of caffeic acid phenylethyl ester in ethanol is approximately 30 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

Caffeic acid phenylethyl ester, an active component of propolis from honeybee hives, is known to have antimutagenic, anticarcinogenic, antiinflammatory, and immunomodulatory properties. It is a potent and specific inhibitor of the activation of nuclear transcription factor NF- κ B.¹ It also exhibits inhibitory activity against HIV-1 integrase with IC₅₀ of 7 μ M.²

References

1. Natarajan, K., Singh, S., Burke, Jr., T.R., *et al.* Caffeic acid phenethyl ester is a potent and specific inhibitor of activation of nuclear transcription factor NF- κ B. *Proc. Natl. Acad. Sci. USA* **93**, 9090-9095 (1996).
2. Burke, Jr., T.R., Fesen, M.R., Mazumder, A., *et al.* Hydroxylated aromatic inhibitors of HIV-1 integrase. *J. Med. Chem.* **38**, 4171-4178 (1995).

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. It is the responsibility of the purchaser to determine suitability for other applications.

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, 08/21/2023

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