

Caffeic acid phenylethyl ester

HIV-1 integrase inhibitor

Active component of propolis from honeybee hives, known to have anti-mitogenic, anti-carcinogenic, anti-inflammatory, and immunomodulatory properties. Potent and specific inhibitor of NF- κ B activation, induced by TNF- α (Prod. No. ALX-522-008 (human) or Prod. No. ALX-522-009 (mouse)), PMA (Prod. No. BML-PE160) and other inflammatory agents. Exhibits inhibitory activity against HIV-1 integrase ($IC_{50}=7\mu M$). Suppresses lipid peroxidation and inhibits ornithine decarboxylase, protein tyrosine kinase and lipoxygenase activities. Induces apoptosis.

Citations: 10

[View Online »](#)

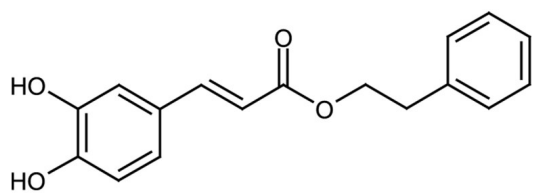
Ordering Information

[Order Online »](#)

ALX-270-244-M010	10mg
ALX-270-244-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored at -20°C.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CAPE, Phenethyl caffeiate, Caffeic acid phenethyl ester
Appearance	White to off-white crystalline solid.
CAS	104594-70-9
Couple Target	HIV
Couple Type	Inhibitor
Formula	$C_{17}H_{16}O_4$
MW	284.3
Purity	≥97% (HPLC)
Solubility	10mg/ml soluble in DMSO, 100% ethanol, methanol, acetone or acetonitrile; insoluble in water.
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)