Caffeic acid phenylethyl ester

HIV-1 integrase inhibitor

Active component of propolis from honeybee hives, known to have antimitogenic, 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 μ M). Suppresses lipid peroxidation and inhibits ornithine decarboxylase, protein tyrosine kinase and lipoxygenase activities. Induces apoptosis.

Citations: 10

View Online »

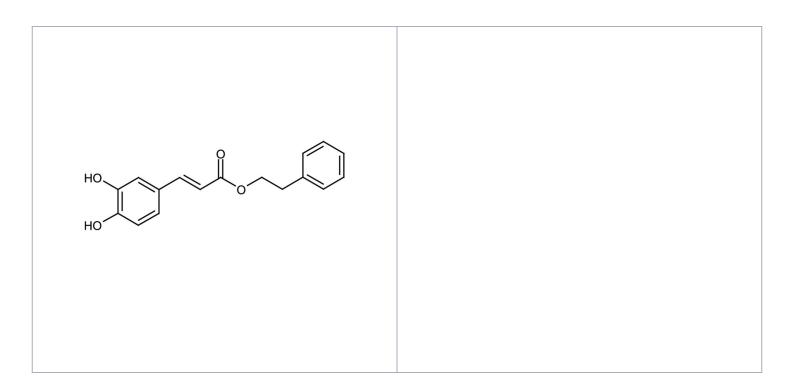
Ordering Information

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 C₁₇H₁₆O₄

MW 284.3

Purity ≥97% (HPLC)

Solubility 10mg/ml soluble in DMSO, 100% ethanol, methanol,

acetone or acetonitrile; insoluble in water.

Source Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com