

8-

Isopentenylaringenin

Phytoestrogen

Prenyl flavonoid. Phytoestrogen. Selective, non-steroidal estrogen receptor α (ER α) ligand. Potent inhibitor of angiogenesis *in vitro* and *in vivo*. Chemopreventive agent against cancer induced by heterocyclic amines.

Citations: 10

[View Online »](#)

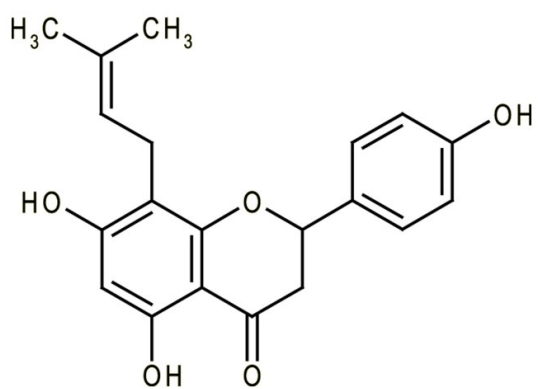
Ordering Information

[Order Online »](#)

ALX-385-025-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Prenylnaringenin

Appearance Light yellow powder

CAS 68682-02-0

Couple Target Estrogen receptor

Couple Type Ligand

Formula $C_{20}H_{20}O_5$

MW 340.4

Purity ≥98% (1H-NMR)

Solubility Soluble in ethanol, methanol or DMSO. Nearly insoluble in water.

Source Semi synthetic, original isolation from hops (*Humulus lupulus* L.)

Last modified: May 29, 2024



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)