

Limonin

Carcinogenesis inhibitor

Bitter principle of citrus fruits. Inhibits chemically induced carcinogenesis. Inhibits HIV-1 replication. Shows antinociceptive effect. Does not possess any inherent antioxidant activity.

Citations: 11

[View Online »](#)

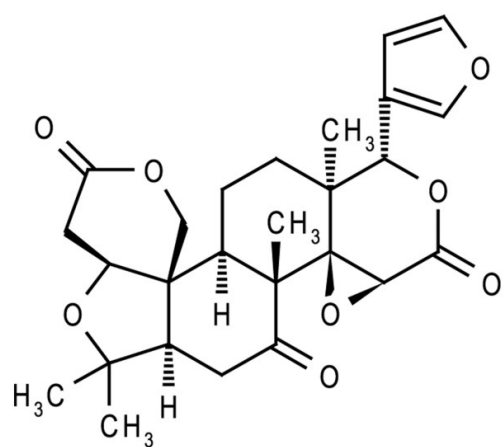
Ordering Information

[Order Online »](#)

ALX-350-225-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Limonoic acid di- δ -lactone

Appearance White to off-white solid.

CAS 1180-71-8

Formula $C_{26}H_{30}O_8$

MI **14:** 5494

MW 470.5

Purity $\geq 98\%$ (UHPLC)

Solubility Soluble in 100% ethanol or glacial acetic acid; slightly soluble in water.

Source Isolated from grapefruit seed.

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-
be@enzolifesciences.com](mailto:info-be@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)