

Xanthohumol

DGAT inhibitor

Prenylated chalcone found in beer. Potent inhibitor of diacylglycerol acetyltransferase (DGAT). Inhibits DNA polymerase and induces cell differentiation. Has antiproliferative and cytotoxic effects in human cancer cell lines. Inhibits human P450 enzymes. Induces quinone reductase. Inhibits the expression of HIF-1 α and VEGF under hypoxic conditions. Induces apoptosis in mature adipocytes.

Citations: 13

[View Online »](#)

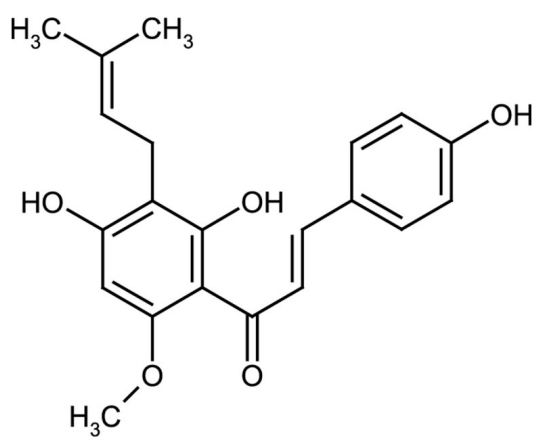
Ordering Information

[Order Online »](#)

ALX-350-280-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solution made in DMSO should not be stored for longer than one month. Working solution is stable for 24 hours at 37°C and up to one week when stored at +4°C. Unstable in acidic (pH 7.5) systems.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow to orange powder.
CAS	569-83-5, 6754-58-1
Couple Target	Cytochrome P450, DGAT, DNA polymerase
Couple Type	Inhibitor
Formula	$C_{21}H_{22}O_5$
Identity	Identity determined by NMR.
MW	354.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), methanol (10mg/ml) or 100% ethanol; soluble in aqueous buffers up to 20μM.
Source	Isolated from hops (<i>Humulus lupulus</i>).

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

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

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