

Ferutinin (high purity)

Estrogen receptor α activator

Potent, naturally occurring non-steroid estrogenic compound. Agonist for estrogen receptor (ER) α and agonist/antagonist for ER β with higher binding affinity than tamoxifen (Prod. No. ALX-550-095) for both ERs. Electrogenic Ca^{2+} ionophore inducing mitochondrial depolarisation which can be completely blocked by cyclosporin A (Prod. No. BML-A195), suggesting that ferutinin opens the mitochondrial permeability transition pore (mPTP). In a concentration range of 1-50 μM ferutinin increases the permeability of thymocytes, mitochondria, sarcoplasmic reticulum, liposomes and bilayer lipid membranes for Ca^{2+} .

Citations: 6

[View Online »](#)

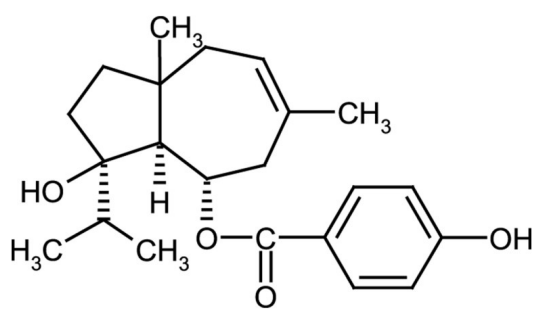
Ordering Information

[Order Online »](#)

ALX-350-098-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

Appearance White to off-white solid

CAS 41743-44-6

Couple Target Estrogen receptor

Couple Type Activator

Formula $C_{22}H_{30}O_4$

MW 358.5

Purity ≥98% (HPLC)

Solubility Soluble in acetone, dichloromethane, DMSO or ethyl acetate.

Source Semisynthetic.

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)