

Orlistat

Fatty acid synthase inhibitor

Cell permeable, irreversible inhibitor of gastric and pancreatic lipases. Shows only minimal activity against amylase, trypsin, chymotrypsin, or phospholipase A₂ (PLA₂). Partially inhibits the hydrolysis of triglycerides and lowers the absorption of dietary fat and promotes weight loss. Anti-obesity drug. Exhibits antitumor activity by inhibition of the thioesterase domain of fatty acid synthase (FAS) both *in vitro* and *in vivo*.

Citations: 8

[View Online »](#)

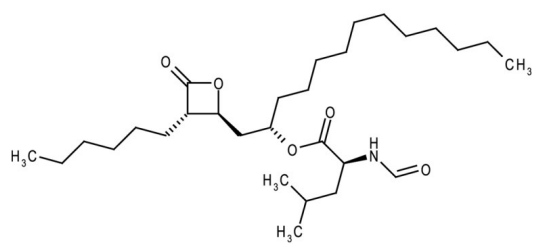
Ordering Information

[Order Online »](#)

ALX-350-152-M250	250mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 6 weeks when stored at -20°C.
Handling	Protect from light and moisture. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tetrahydrolipstatin, N-Formyl-L-leucine-(1S)-1-(((2S,3S)-3-hexyl-4-oxo-2-oxetanyl)methyl)dodecyl ester
Appearance	White to off-white solid.
CAS	96829-58-2
Couple Target	Fatty acid synthase
Couple Type	Inhibitor
Formula	$C_{29}H_{53}NO_5$
MI	14: 6869
MW	495.7
Purity	≥97%
Solubility	Soluble in DMSO or 100% ethanol.
Source	Synthetic. Originally isolated from <i>Streptomyces</i> sp.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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