

C75

Fatty acid synthase inhibitor

Inhibitor of fatty acid synthase (FAS) reducing food intake and body weight in mice. Exhibits irreversible slow-binding biphasic inactivation of fatty acid synthase. Downregulates neuropeptide Y and Agouti-related protein expression. Has been proposed to activate CPT-1 activity in liver and adipose tissue, leading to increased fatty acid oxidation and energy production. Shows significant *in vivo* antitumor activity in human breast cancer cells. Suppresses DNA replication and induces apoptosis. FAS inhibition by C75 leads to a dramatic accumulation of the CDK inhibitor p27KIP1 from cytosol to cell nuclei.

Citations: 26

[View Online »](#)

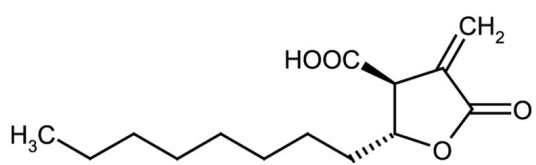
Ordering Information

[Order Online »](#)

ALX-270-286-M001	1mg
ALX-270-286-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	trans-4-Carboxy-5-octyl-3-methylene-butyrolactone
Appearance	White to off-white solid.
CAS	191282-48-1
Couple Target	Fatty acid synthase
Couple Type	Inhibitor
Formula	$C_{14}H_{22}O_4$
MW	254.3
Purity	≥98%
Solubility	Soluble in dichloromethane, methanol or DMSO.
Source	Synthetic.



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)