

Amidepsine A

DGAT inhibitor

Inhibitor of diacylglycerol acyltransferase (DGAT). Shows inhibitory effect in rat liver microsomes ($IC_{50}=10.2\mu\text{M}$) and Raji cells ($IC_{50}=15.5\mu\text{M}$). Excessive accumulation of triacetylglycerol produced by DGAT may cause fatty liver, obesity and hypertriglyceridemia, which may lead to atherosclerosis, diabetes and metabolic disorders.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-380-262-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Pale white to pale yellow powder.

CAS 169181-28-6

Couple Target DGAT

Couple Type Inhibitor

Formula $C_{29}H_{29}NO_{11}$

MW 567.6

Purity ≥95% (HPLC)

Solubility Soluble in methanol, ethyl acetate or benzene; insoluble in water or hexane.

Source Isolated from fungal strain FO-2943.

Last modified: May 29, 2024



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-be@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