

Clofibrate

PPAR α activator

Activates PPAR α . Induces cytochrome P450 4A1 and 4A3 which catalize the omega-oxidation of fatty acids and prostaglandins.

Citations: 3

[View Online »](#)

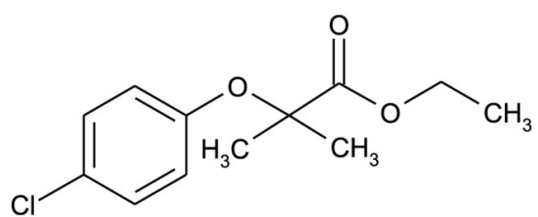
Ordering Information

[Order Online »](#)

ALX-270-256-M500	500mg
ALX-270-256-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for up to 6 months when stored at –20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(p-Chlorophenoxy)-2-methylpropionic acid ethyl ester
Appearance	Colorless oil.
CAS	637-07-0
Couple Target	PPAR
Couple Type	Activator, Ligand
Formula	$C_{12}H_{15}ClO_3$
MI	14: 2377
MW	242.7
Purity	≥98% (GC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)