

Ciprofibrate

PPAR α activator

Hypolipidemic compound. PPAR α (peroxisome proliferator-activated receptor α) agonist.

Citations: 4

[View Online »](#)

Ordering Information

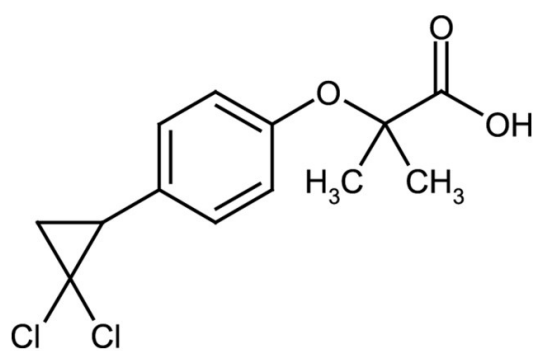
[Order Online »](#)

ALX-270-475-M025

25mg

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

Alternative Name 2-[4-(2,2-Dichlorocyclopropyl)phenoxy]-2-methylpropanoic acid

Appearance White to light yellow powder.

CAS 52214-84-3

Couple Target PPAR

Couple Type Activator, Ligand

Formula $C_{13}H_{14}Cl_2O_3$

MI **14:** 2313

MW 289.2

Purity ≥98% (HPLC)

RTECS UF0880000

Solubility Soluble in DMSO.

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-
fr@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)