

Fenofibrate

PPAR α activator

Hypolipidemic compound. PPAR α (peroxisome proliferator-activated receptor α) agonist (EC_{50} =18 μ M for mouse and EC_{50} =30 μ M for human PPAR α). Also binds to PPAR γ , but with at least 10-fold less affinity and is inactive at PPAR δ (up to EC_{50} =100 μ M).

Citations: 6

[View Online »](#)

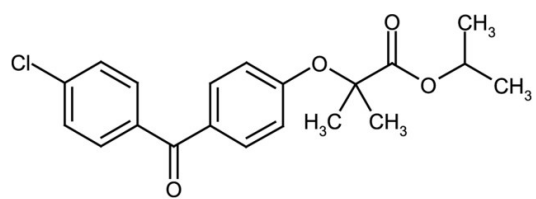
Ordering Information

[Order Online »](#)

ALX-270-481-G005	5g
------------------	----

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-(4-Chlorobenzoyl)phenoxy]-2-methyl-propanoic acid
1-methylethyl ester, Procetofen

Appearance White to off-white crystalline solid.

CAS 49562-28-9

Couple Target PPAR

Couple Type Activator, Ligand

Formula $C_{20}H_{21}ClO_4$

MI 14: 3978

MW 360.8

Purity ≥98% (GC)

RTECS UA2453400

Solubility Soluble in acetone, ether, DMSO or dimethyl formamide; slightly soluble in 100% ethanol or methanol; insoluble in water.

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)