

WY-14643

PPAR ligand

WY-14643 is a potent and widely used peroxisome proliferator activator receptor (PPAR) activator. It has been shown to be selective for PPAR α .

Citations: 14

[View Online »](#)

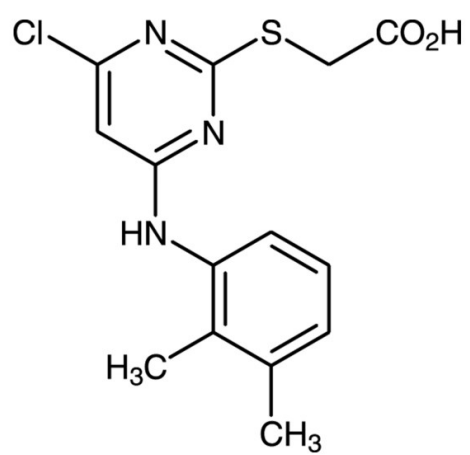
Ordering Information

[Order Online »](#)

BML-GR200-0250	250mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Pirinixic acid

Appearance White to off-white solid.

CAS 50892-23-4

Couple Target PPAR

Couple Type Ligand

Formula $C_{14}H_{14}ClN_3O_2S$

Identity Determined by NMR.

MW 323.8

MeltingPoint 155-157°C

Purity ≥97% (HPLC)

RTECS AG2915000

Solubility Soluble in ethyl acetate or dilute base or slightly soluble in alcohol.

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)