

Rosiglitazone

Insulin sensitizing agent

Potent insulin sensitizing agent. Peroxisome proliferator-activated receptor γ (PPAR- γ) agonist. Hepatotoxic. Antidiabetic.

Citations: 51

[View Online »](#)

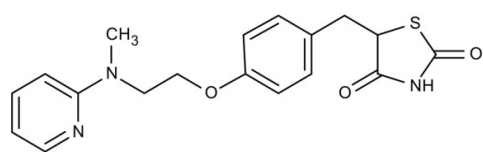
Ordering Information

[Order Online »](#)

ALX-350-125-M025	25mg
ALX-350-125-M100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 2 years after receipt when stored at +4°C. Stock solutions are stable for up to 3 months when stored at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	122320-73-4
Couple Target	PPAR
Couple Type	Activator, Ligand
Formula	$C_{18}H_{19}N_3O_3S$
MI	14: 8265
MW	357.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml), 100% ethanol or methanol.

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)