

Pioglitazone

PPAR γ activator

Pioglitazone selectively activates PPAR γ -1. It is about one tenth as potent as rosiglitazone (EC_{50} ~500nM for human and mouse PPAR γ). In a transactivation assay using COS-1 cells transfected with full length human PPAR α and RXR α , pioglitazone and rosiglitazone exhibit low level activation of PPAR α at 1 μ M and 5.4- and 4.2-fold activation, respectively, at a concentration of 10 μ M.

Thiazolidinediones (TZDs) are a group of structurally related PPAR γ agonists with anti-diabetic actions in vivo. Rosiglitazone (Prod. No. ALX-350-125) is a prototypical TZD and has served as a reference compound for this class of PPAR γ ligands.

Citations: 13

[View Online »](#)

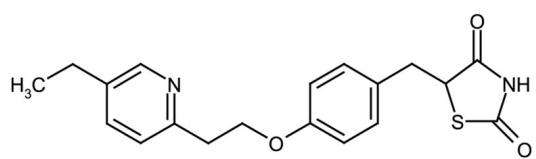
Ordering Information

[Order Online »](#)

ALX-270-367-M001	1mg
ALX-270-367-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	111025-46-8
Couple Target	PPAR
Couple Type	Activator, Ligand
Formula	$C_{19}H_{20}N_2O_3S$
Identity	Confirmed by 1H-NMR
MI	14: 7452
MW	356.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (2.5mg/ml) or dimethyl formamide (2.5mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com