

Ciglitazone

PPAR ligand

Ciglitazone is a thiazolidinedione type of antihyperglycemic agent. It displays antihyperglycemic activity in genetically obese C57 Bl/6 *ob/ob* mice and is a selective PPAR γ agonist (EC_{50} =3 μ M). Inhibits HUVEC differentiation and angiogenesis and also stimulates adipogenesis and decreases osteoblastogenesis in human mesenchymal stem cells.

Citations: 8

[View Online »](#)

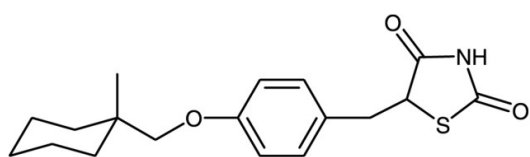
Ordering Information

[Order Online »](#)

BML-GR205-0005	5mg
BML-GR205-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white.
CAS	74772-77-3
Couple Target	PPAR
Couple Type	Ligand
Formula	$C_{18}H_{23}NO_3S$
Identity	Determined by NMR.
MW	333.5
Purity	≥98% (HPLC)
RTECS	XJ5813700
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).
Source	Synthetic.

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)