

Troglitazone

PPAR ligand

Antidiabetic thiazolidinedione. Potent and selective PPAR γ ligand. Binds to the PPAR γ ligand-binding domain and recruits ligand-specific transcriptional coactivators at 1 μ M. In skeletal muscle and adipocyte cell cultures, the EC₅₀ of troglitazone is about 10 μ M. Displaces [125I]SB-236636 with an IC₅₀=7.9 μ M. Stimulates adipogenesis and decreases osteoblastogenesis in human mesenchymal stem cells. Induces apoptosis in human vascular smooth muscle cells and inhibits the growth of a variety of human cancer cells including lung, bladder and myeloid leukemia cells. Displays an anti-invasive effect on human breast cancer cells; reduces migration, adhesion and spreading on fibronectin-coated plates. Inhibits cell growth of hematopoietic cell lines. Inhibits lamellipodia formation and actin polymerization.

Citations: 12

[View Online »](#)

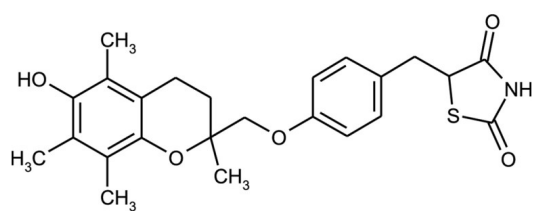
Ordering Information

[Order Online »](#)

BML-GR210-0005	5mg
BML-GR210-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-5-[[4-[(3,4-Dihydro-6-hydroxy-2,5,7,8-tetramethyl-2H-1-benzopyran-2-yl)methoxy]phenyl]methyl]-2,4-thiazolidinedione
Appearance	White to yellow solid.
CAS	97322-87-7
Couple Target	PPAR
Couple Type	Ligand
Formula	C ₂₄ H ₂₇ NO ₅ S
MI	14: 9769
MW	441.5
Purity	≥98% (HPLC)
RTECS	XJ5813130
Solubility	Soluble in DMSO (>25mg/ml), 100% ethanol (>25mg/ml, warm) or dimethyl formamide.
Source	Synthetic.

Technical Info / Product Notes

Troglitazone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, troglitazone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Troglitazone has a solubility of 100µg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Note: Product is not sterile.



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-
be@enzolifesciences.com](mailto:info-be@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)