# **Troglitazone**

**PPAR ligand** 

Antidiabetic thiazolidinedione. Potent and selective PPAR $\gamma$  ligand. Binds to the PPAR $\gamma$  ligand-binding domain and recruits ligand-specific transcriptional coactivators at 1 $\mu$ M. In skeletal muscle and adipocyte cell cultures, the EC $_{50}$  of troglitazone is about 10 $\mu$ M. Displaces [125I]SB-236636 with an IC $_{50}$ =7.9 $\mu$ 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 »

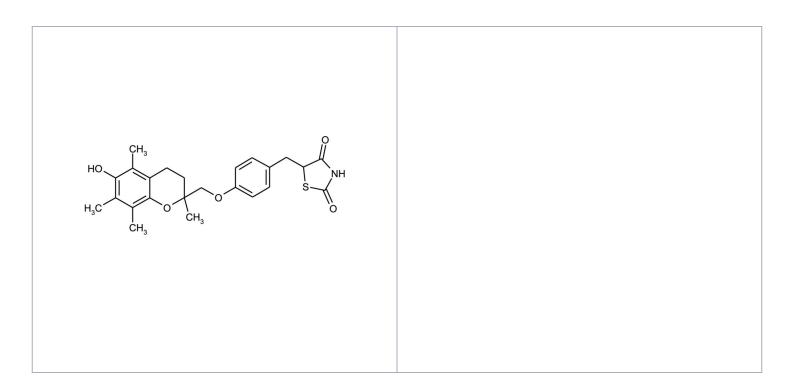
## **Ordering Information**

**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  $C_{24}H_{27}NO_5S$ 

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.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com