

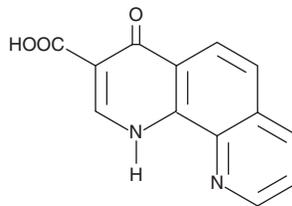
# PRODUCT INFORMATION



## Ciglitazone

Item No. 71730

**CAS Registry No.:** 74772-77-3  
**Formal Name:** 5-[[4-[(1-methylcyclohexyl) methoxy]phenyl]methyl]-2,4-thiazolidinedione  
**Synonyms:** ADD 3878, U-63287  
**MF:** C<sub>18</sub>H<sub>23</sub>NO<sub>3</sub>S  
**FW:** 333.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 228, 279, 284 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Ciglitazone is supplied as a crystalline solid. A stock solution may be made by dissolving the ciglitazone in the solvent of choice, which should be purged with an inert gas. Ciglitazone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ciglitazone in these solvents is approximately 29, 21, and 16 mg/ml, respectively.

Ciglitazone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Ciglitazone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ciglitazone has a solubility of approximately 400 µg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Ciglitazone is an antidiabetic drug of the thiazolidinedione structural class. Ciglitazone is a potent and selective PPAR $\gamma$  ligand. It binds to the PPAR $\gamma$  ligand-binding domain with an EC<sub>50</sub> of 3.0 µM.<sup>1</sup> Ciglitazone is active *in vivo* as an anti-hyperglycemic agent in the *ob/ob* mouse model.<sup>1</sup>

### Reference

1. Willson, T.M., Cobb, J.E., Cowan, D.J., *et al.* The structure-activity relationship between peroxisome proliferator-activated receptor  $\gamma$  agonism and the antihyperglycemic activity of thiazolidinediones. *J. Med. Chem.* **39**(3), 665-668 (1996).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/08/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM