

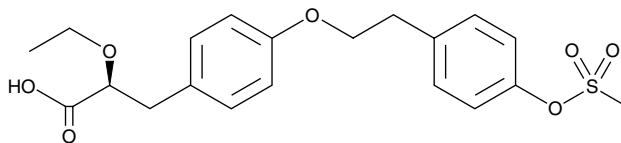
# Product Information



## Tesaglitazar

Item No. 16791

**CAS Registry No.:** 251565-85-2  
**Formal Name:** ( $\alpha$ S)-ethoxy-4-[2-[4-[(methylsulfonyl)oxy]phenyl]ethoxy]-benzenepropanoic acid  
**Synonyms:** AZ 242, Galida™  
**MF:** C<sub>20</sub>H<sub>24</sub>O<sub>7</sub>S  
**FW:** 408.5  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 225, 275 nm



### Laboratory Procedures

For long term storage, we suggest that tesaglitazar be stored as supplied at -20°C. It should be stable for at least two years.

Tesaglitazar is supplied as a crystalline solid. A stock solution may be made by dissolving the tesaglitazar in the solvent of choice. Tesaglitazar is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of tesaglitazar in ethanol is approximately 10 mg/ml and approximately 50 mg/ml in DMSO and DMF.

Tesaglitazar is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tesaglitazar should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Tesaglitazar has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Peroxisome proliferator-activated receptors (PPARs) are activated by fatty acids and eicosanoids as well as antidyslipidemic agents. Among the receptor isotypes, PPAR $\alpha$  demonstrates a particular role in fatty acid oxidation whereas PPAR $\gamma$  is known to be involved in adipocyte differentiation and lipid storage. Tesaglitazar, a dihydro cinnamate derivative, is a dual agonist of PPAR $\alpha$  and  $\gamma$  that demonstrates IC<sub>50</sub> values of 1 and 0.2  $\mu$ M, respectively in ligand binding assays.<sup>1</sup> At 3  $\mu$ M/kg/day for three weeks, tesaglitazar has been used to reduce insulin resistance in obese Zucker rats.<sup>2</sup> Furthermore, it has been investigated clinically for its potential to address disorders in glucose and lipid metabolism in patients with type 2 diabetes.<sup>3</sup>

### References

1. Cronet, P., Petersen, J.F.W., Folmer, R., *et al.* Structure of the PPAR $\alpha$  and - $\gamma$  ligand binding domain in complex with AZ 242; ligand selectivity and agonist activation in the PPAR family. *Structure* **9(8)**, 699-706 (2001).
2. Wallenius, K., Kjellstedt, A., Thalén, P., *et al.* The PPAR $\alpha$ / $\gamma$  agonist, tesaglitazar, improves insulin mediated switching of tissue glucose and free fatty acid utilization *in vivo* in the obese Zucker rat. *PPAR Res.* **2013**, 1-14 (2013).
3. Wilding, J.P.H., Gause-Nilsson, I., Persson, A., *et al.* Tesaglitazar, as add-on therapy to sulphonylurea, dose-dependently improves glucose and lipid abnormalities in patients with type 2 diabetes. *Diab. Vasc. Dis. Res.* **4(3)**, 194-203 (2007).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/16791](http://www.caymanchem.com/catalog/16791)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 11/05/2014

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)