

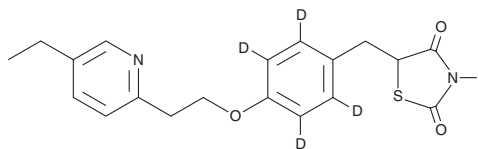
# PRODUCT INFORMATION



## Pioglitazone-d<sub>4</sub>

Item No. 18259

**CAS Registry No.:** 1134163-29-3  
**Formal Name:** 5-[[4-[2-(5-ethyl-2-pyridinyl)ethoxy]phenyl-2,3,5,6-d<sub>4</sub>]methyl]-2,4-thiazolidinedione  
**MF:** C<sub>19</sub>H<sub>16</sub>D<sub>4</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 360.5  
**Chemical Purity:** ≥98% (Pioglitazone)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A 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

Pioglitazone-d<sub>4</sub> is intended for use as an internal standard for the quantification of pioglitazone (Item Nos. 71745 | 10028 | 22263) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Pioglitazone-d<sub>4</sub> is supplied as a solid. A stock solution may be made by dissolving the pioglitazone-d<sub>4</sub> in the solvent of choice. Pioglitazone-d<sub>4</sub> is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of pioglitazone-d<sub>4</sub> in these solvents is approximately 2.5 mg/ml.

### Description

Pioglitazone is an agonist of the peroxisome proliferator-activated receptor  $\gamma$  (PPAR $\gamma$ ; EC<sub>50</sub> = ~500-600 nM for both human and murine PPAR $\gamma$ ).<sup>1,2</sup> It is selective for PPAR $\gamma$  over PPAR $\alpha$ , exhibiting low level activation of PPAR $\alpha$  at 1  $\mu$ M and 5.4-fold activation at a concentration of 10  $\mu$ M.<sup>1</sup> Pioglitazone inhibits pyruvate oxidation and glucose production in hepatocytes when used at a concentration of 10  $\mu$ M.<sup>3</sup> *In vivo*, pioglitazone (0.3-3 mg/kg per day) reduces hyperglycemia, hyperlipidemia, and hyperinsulinemia in a dose-dependent manner in male Wistar fatty rats.<sup>4</sup> It reduces the number of lesions in a transgenic rat adenocarcinoma of prostate (TRAP) model.<sup>5</sup> Pioglitazone (2.5 mg/kg) also decreases production of neuroinflammatory cytokines and reduces immobility in the forced swim and tail suspension tests in a mouse model of chronic mild stress, indicating antidepressant-like activity that can be reversed by the PPAR $\gamma$  antagonist GW 9662 (Item No. 70785).<sup>6</sup>

### References

1. Sakamoto, J., Kimura, H., Moriyama, S., et al. *Biochem. Biophys. Res. Commun.* **278**(3), 704-711 (2000).
2. Willson, T.M., Brown, P.J., Sternbach, D.D., et al. *J. Med. Chem.* **43**(4), 528-550 (2000).
3. Shannon, C.E., Daniele, G., Galindo, C., et al. *FEBS J.* **284**(3), 451-465 (2017).
4. Sugiyama, Y., Taketomi, S., Shimura, Y., et al. *Arzneimittelforschung.* **40**(3), 263-267 (1990).
5. Suzuki, S., Mori, Y., Nagano, A., et al. *Int. J. Mol. Sci.* **17**(12), pii: E2071 (2016).
6. Zhao, Q., Wu, X., Yan, S., et al. *J. Neuroinflammation* **13**(1), 259 (2016).

#### 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, 12/08/2022

#### 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