

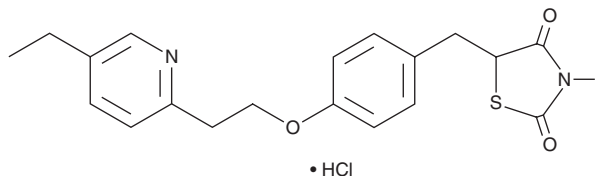
PRODUCT INFORMATION



Pioglitazone (hydrochloride)

Item No. 22263

CAS Registry No.: 112529-15-4
Formal Name: 5-[[4-[2-(5-ethyl-2-pyridinyl)ethoxy]phenyl]methyl]-2,4-thiazolidinedione, monohydrochloride
MF: C₁₉H₂₀N₂O₃S • HCl
FW: 392.9
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 269 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

Pioglitazone (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the pioglitazone (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Pioglitazone (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pioglitazone (hydrochloride) in these solvents is approximately 20 mg/ml.

Pioglitazone (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pioglitazone (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pioglitazone (hydrochloride) has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pioglitazone is an agonist of the peroxisome proliferator-activated receptor γ (PPAR γ ; EC₅₀ = ~500-600 nM for both human and murine PPAR γ).^{1,2} It is selective for PPAR γ over PPAR α , exhibiting low level activation of PPAR α at 1 μ M and 5.4-fold activation at a concentration of 10 μ M.¹ Pioglitazone inhibits pyruvate oxidation and glucose production in hepatocytes when used at a concentration of 10 μ M.³ *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.⁴ It reduces the number of lesions in a transgenic rat adenocarcinoma of prostate (TRAP) model.⁵ 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 γ antagonist GW 9662 (Item No. 70785).⁶

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), 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, 12/19/2022

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