

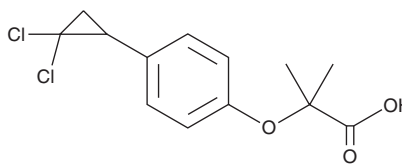
PRODUCT INFORMATION



Ciprofibrate

Item No. 18515

CAS Registry No.: 52214-84-3
Formal Name: 2-[4-(2,2-dichlorocyclopropyl)phenoxy]-2-methyl-propanoic acid
Synonyms: (±)-Ciprofibrate, WIN 35,833
MF: C₁₃H₁₄Cl₂O₃
FW: 289.2
Purity: ≥98%
UV/Vis.: λ_{max}: 232 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

Ciprofibrate is supplied as a crystalline solid. A stock solution may be made by dissolving the ciprofibrate in the solvent of choice, which should be purged with an inert gas. Ciprofibrate is soluble in the organic solvent methanol at a concentration of approximately 20 mg/ml.

Description

Ciprofibrate is an agonist of peroxisome proliferator-activated receptor α (PPAR α ; EC₅₀ = 0.9 μ M in a transactivation assay).¹ It is selective for PPAR α over PPAR γ and PPAR δ at 300 μ M.² Ciprofibrate (250 μ M) induces cell cycle arrest at the G₂/M and S phases in Fao rat, but not HepG2 human, hepatocellular carcinoma cells.³ It decreases fasting plasma levels of triglycerides and increases fasting plasma glucose levels in the apolipoprotein CIII transgenic mouse model of hypertriglyceridemia when administered at a dose of 10 mg/kg.⁴ Formulations containing ciprofibrate have been used in the treatment of hypertriglyceridemia.

References

1. Quang, T.H., Ngan, N.T.T., Minh, C.V., *et al.* Anti-inflammatory and PPAR transactivational effects of secondary metabolites from the roots of *Asarum sieboldii*. *Bioorg. Med. Chem. Lett.* **22(7)**, 2527-2533 (2012).
2. Forman, B.M., Chen, J., and Evans, R.M. Hypolipidemic drugs, polyunsaturated fatty acids, and eicosanoids are ligands for peroxisome proliferator-activated receptors α and δ . *Proc. Natl. Acad. Sci. USA* **94(9)**, 4312-4317 (1997).
3. Passilly, P., Jannin, B., Hassell, S.J., *et al.* Human HepG2 and rat Fao hepatic-derived cell lines show different responses to ciprofibrate, a peroxisome proliferator: analysis by flow cytometry. *Exp. Cell. Res.* **223(2)**, 436-442 (1996).
4. Bighetti, E.J.B., Patricio, P.R., Casquero, A.C., *et al.* Ciprofibrate increases cholesteryl ester transfer protein gene expression and the indirect reverse cholesterol transport to the liver. *Lipids Health Dis.* **8**, 50 (2009).

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, 01/03/2023

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