

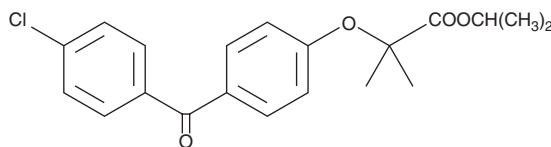
# PRODUCT INFORMATION



## Fenofibrate

Item No. 10005368

**CAS Registry No.:** 49562-28-9  
**Formal Name:** 2-[4-(4-chlorobenzoyl)phenoxy]-2-methyl-propanoic acid, 1-methylethyl ester  
**MF:** C<sub>20</sub>H<sub>21</sub>ClO<sub>4</sub>  
**FW:** 360.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 287 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

Fenofibrate is supplied as a crystalline solid. A stock solution may be made by dissolving the fenofibrate in the solvent of choice, which should be purged with an inert gas. Fenofibrate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fenofibrate in these solvents is 1, 15, and 30 mg/ml, respectively.

Fenofibrate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fenofibrate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Fenofibrate has a solubility of 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Fenofibrate is an agonist of peroxisome proliferator-activated receptor  $\alpha$  (PPAR $\alpha$ ) with EC<sub>50</sub> values of 18 and 30  $\mu$ M for mouse and human receptors, respectively, in a transactivation assay.<sup>1</sup> It is selective for PPAR $\alpha$  over PPAR $\gamma$  (EC<sub>50</sub>s = 300 and 200  $\mu$ M for mouse and human receptors, respectively) and lacks activity at mouse and human PPAR $\delta$  at a concentration of 100  $\mu$ M. *In vivo*, fenofibrate (50-100 mg/kg) reduces plasma levels of triglycerides, C-reactive protein, and malondialdehyde (MDA) in mice with fructose-induced hypertriglyceremia in a dose-dependent manner.<sup>2</sup> It decreases glomerular and tubular atrophy and necrosis induced by cisplatin (Item No. 13119) in rat kidney when administered at a dose of 100 mg/kg.<sup>3</sup> Fenofibrate also reduces the number of pulmonary lesions induced by 4-nitroquinoline 1-oxide (4-NQO) in lung in Tsumura Suzuki obese diabetic (TSOD) mice.<sup>4</sup>

### References

- Willson, T.M., Brown, P.J., Sternbach, D.D., *et al.* The PPARs: From orphan receptors to drug discovery. *J. Med. Chem.* **43**(4), 528-550 (2000).
- Sun, B., Xie, Y., Jiang, J., *et al.* Pleiotropic effects of fenofibrate therapy on rats with hypertriglyceremia. *Lipids Health Dis.* **14**:27, (2015).
- Helmy, M.M., Helmy, M.W., and El-Mas, M.M. Additive renoprotection by pioglitazone and fenofibrate against inflammatory, oxidative and apoptotic manifestations of cisplatin nephrotoxicity: Modulation by PPARs. *PLoS One* **10**(11), e0142303 (2015).
- Kuno, T., Hata, K., Takamatsu, M., *et al.* The peroxisome proliferator-activated receptor (PPAR)  $\alpha$  agonist fenofibrate suppresses chemically induced lung alveolar proliferative lesions in male obese hyperlipidemic mice. *Int. J. Mol. Sci.* **15**(5), 9160-9172 (2014).

#### 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, 11/10/2022

#### 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