

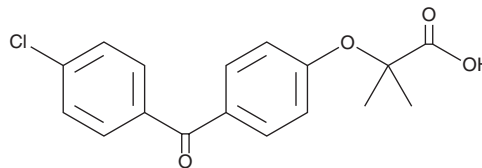
# PRODUCT INFORMATION



## Fenofibric Acid

Item No. 19262

**CAS Registry No.:** 42017-89-0  
**Formal Name:** 2-[4-(4-chlorobenzoyl)phenoxy]-2-methylpropanoic acid  
**Synonyms:** FNF Acid, NSC 281318, Procetofenic Acid  
**MF:** C<sub>17</sub>H<sub>15</sub>ClO<sub>4</sub>  
**FW:** 318.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 288 nm  
**Supplied as:** A crystalline solid  
**Storage:** 4°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Fenofibric acid is supplied as a crystalline solid. A stock solution may be made by dissolving the fenofibric acid in the solvent of choice. Fenofibric acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of fenofibric acid in these solvents is approximately 16, 2, and 14 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of fenofibric acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of fenofibric acid in PBS, pH 7.2, is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Fenofibric acid is an agonist of peroxisome proliferator-activated receptor  $\alpha$  (PPAR $\alpha$ ), PPAR $\gamma$ , and PPAR $\delta$  (EC<sub>50</sub>s = 22.4, 1.47, and 1.06  $\mu$ M, respectively, in a transactivation assay) and an active metabolite of the PPAR $\alpha$  agonist fenofibrate (Item No. 10005368).<sup>1,2</sup> It is formed by the hydrolysis of fenofibrate in tissue.<sup>1</sup> It inhibits COX-2 activity in a cell-free assay (IC<sub>50</sub> = 0.048  $\mu$ M).<sup>4</sup> Fenofibric acid (500  $\mu$ M) reduces aldehyde oxidase 1 (AOX1) protein levels in HepG2 cells.<sup>3</sup> Fenofibric acid (10 mg/kg) reduces carrageenan-induced paw edema in rats.<sup>4</sup> Formulations containing fenofibric acid have been used in the treatment of high cholesterol and hypertriglyceridemia.

### References

1. Caldwell, J. The biochemical pharmacology of fenofibrate. *Cardiology* **76(Suppl 1)**, 33-44 (1989).
2. Dietz, M., Mohr, P., Kuhn, B., et al. Comparative molecular profiling of the PPAR $\alpha$ / $\gamma$  activator aleglitazar: PPAR selectivity, activity and interaction with cofactors. *ChemMedChem* **7(6)**, 1101-1111 (2012).
3. Neumeier, M., Weigert, J., Schäffler, A., et al. Aldehyde oxidase 1 is highly abundant in hepatic steatosis and is downregulated by adiponectin and fenofibric acid in hepatocytes *in vitro*. *Biochem. Biophys. Res. Commun.* **350(3)**, 731-735 (2006).
4. Prasad, G.S., Govardhan, P., Deepika, G., et al. Anti-inflammatory activity of anti-hyperlipidemic drug, fenofibrate, and its phase-I metabolite fenofibric acid: *in silico*, *in vitro*, and *in vivo* studies. *Inflammopharmacology* **26(4)**, 973-981 (2018).

#### 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, 11/11/2022

#### 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