

# PRODUCT INFORMATION



**CP 775,146**

Item No. 28290

**CAS Registry No.:** 702680-17-9

**Formal Name:** 2-methyl-2-[3-[(3S)-1-[2-[4-(1-methylethyl)phenyl]acetyl]-3-piperidinyl]phenoxy]-propanoic acid

**MF:** C<sub>26</sub>H<sub>33</sub>NO<sub>4</sub>

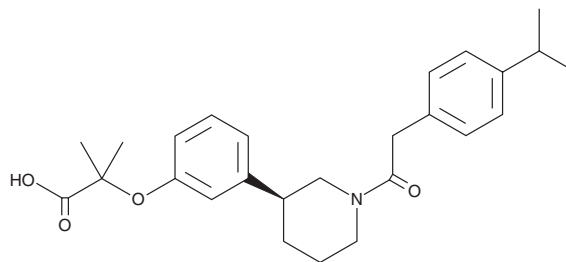
**FW:** 423.5

**Purity:** ≥95%

**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

CP 775,146 is supplied as a solid. A stock solution may be made by dissolving the CP 775,146 in the solvent of choice, which should be purged with an inert gas. CP 775,146 is slightly soluble in methanol.

## Description

CP 775,146 is an agonist of peroxisome proliferator-activated receptor  $\alpha$  (PPAR $\alpha$ ) that selectively binds to PPAR $\alpha$  ( $K_i = 24.5$  nM) over PPAR $\beta$  and PPAR $\gamma$  ( $K_{iS} = >10,000$  nM).<sup>1</sup> CP 775,146 induces PPAR $\alpha$  transcriptional activity in a reporter assay ( $EC_{50S} = 57$  and 284 nM for human and rat receptors, respectively, expressed in HepG2 cells). It decreases plasma triglyceride levels in mice by 73% when administered at a dose of 2 mg/kg per day for two days.

## Reference

1. Kane, C.D., Stevens, K.A., Fischer, J.E., *et al.* Molecular characterization of novel and selective peroxisome proliferator-activated receptor alpha agonists with robust hypolipidemic activity in vivo. *Mol. Pharmacol.* 75(2), 296-306 (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, 10/26/2022

## 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