# **PRODUCT** INFORMATION



GW 590735

Item No. 10009880

CAS Registry No.:	343321-96-0	
Formal Name:	2-methyl-2-[4-[[[[4-methyl-2-[4-	
	(trifluoromethyl)phenyl]-5-thiazolyl]carbonyl]	
	amino]methyl]phenoxy]-propanoic acid	0
MF:	$C_{23}H_{21}F_{3}N_{2}O_{4}S$	
FW:	478.5	
Purity:	≥ <b>98%</b> н	
UV/Vis.:	λ <sub>max</sub> : 224, 311 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	
Stability:	≥2 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis		

# Laboratory Procedures

GW 590735 is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 590735 in the solvent of choice, which should be purged with an inert gas. GW 590735 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GW 590735 in these solvents is approximately 30 mg/ml.

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 GW 590735 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of GW 590735 in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

# Description

The peroxisome proliferator-activated receptors (PPARs)  $\alpha$ ,  $\gamma$ , and  $\delta$  are ligand-activated transcription factors that play a key role in lipid homeostasis. Activation of PPARa results in increased clearance of triglyceride (TG) rich very low-density lipoprotein (VLDL) via a reduction in plasma levels of ApoCIII and in upregulation of ApoA1, the principal lipoprotein component of high-density lipoprotein (HDL).<sup>1</sup> GW 590735 is a potent and selective agonist of PPAR $\alpha$  with an EC<sub>50</sub> value of 4 nM for the expression of a GAL4-responsive reporter gene and at least 500-fold selectivity versus PPARy and PPARo.<sup>1</sup>

# Reference

1. Sierra, M.L., Beneton, V., Boullay, A.-B., et al. Subsituted 2-[(4-aminomethyl)phenoxy]-2-methylpropionic acid PPARα agonists. 1. Discovery of a novel series of potent HDLc raising agents. J. Med. Chem. 50(4), 685-695 (2007).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

## SAFFTY 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, 11/15/2023

# 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