

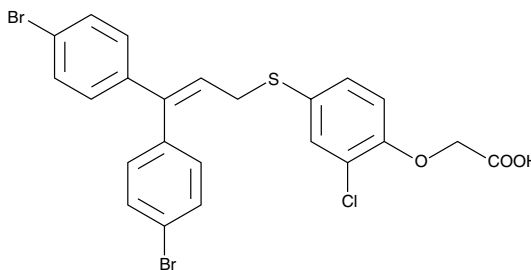
Product Information



CAY10592

Item No. 10012536

CAS Registry No.: 685139-10-0
Formal Name: 2-[4-[[3,3-bis(4-bromophenyl)-2-propen-1-yl]thio]-2-chlorophenoxy]-acetic acid
MF: C₂₃H₁₇Br₂ClO₃S
FW: 568.7
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 243 nm



Laboratory Procedures

For long term storage, we suggest that CAY10592 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10592 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10592 in an organic solvent purged with an inert gas. CAY10592 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10592 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CAY10592 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10592 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10592 has a solubility of approximately 0.2 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Peroxisome proliferator-activated receptors (PPARs) α , δ , γ are ligand-activated nuclear transcription factors involved in the regulation of energy homeostasis as well as insulin sensitivity and glucose metabolism.^{1,2} Pharmacologies of PPAR δ receptor agonists, though relatively obscure, have recently been reported to elevate high-density lipoprotein (HDL) cholesterol and lower plasma triglyceride (TG) levels in obese insulin resistant rhesus monkeys.³ CAY10592 is a full PPAR δ agonist (EC₅₀ = 30 nM) in a fatty acid oxidation assay of rat L6 muscle cells with desirable oral pharmacokinetic properties.⁴ In a transactivation assay using human PPAR receptors, CAY10592 acts as a selective partial PPAR δ agonist (EC₅₀ = 53 nM) with no effect on PPAR α or PPAR γ activity up to 30 μ M.⁴ Chronic treatment of high fat fed ApoB100/CETP-transgenic mice with CAY10592 at a dose of 20 mg/kg increases HDL levels, decreases low-density lipoprotein and TG levels, and improves insulin sensitivity.⁴

References

1. Michalik, L., Auwerx, J., Berger, J.P., *et al.* International union of pharmacology. LXI. Peroxisome proliferator-activated receptors. *Pharmacol. Rev.* **58**, 726-741 (2006).
2. Bratton, L.D., Filzen, G.F., Geyer, A., *et al.* Discovery of highly potent and selective benzyloxybenzyl-based peroxisome proliferator-activator receptor (PPAR) δ agonists. *Bioorganic & Medicinal Chemistry Letters* **17**, 3624-3629 (2007).
3. Oliver, W.R., Shenk, J.L., Snaith, M.R., *et al.* A selective peroxisome proliferator-activated receptor δ agonist promotes reverse cholesterol transport. *Proc. Natl. Acad. Sci. USA* **98(9)**, 5306-5311 (2001).
4. Sauerberg, P., Olsen, G.S., Jeppesen, L., *et al.* Identification and synthesis of a novel selective partial PPAR δ agonist with full efficacy on lipid metabolism *in vitro* and *in vivo*. *J. Med. Chem.* **50**, 1495-1503 (2007).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10012536

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 05/02/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com