

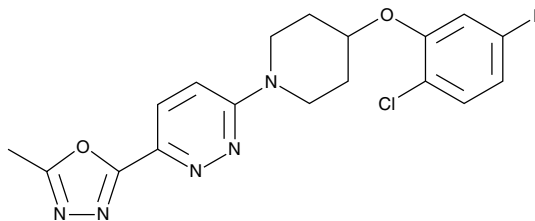
Product Information



CAY10566

Item No. 10012562

CAS Registry No.: 944808-88-2
Formal Name: 3-[4-(2-chloro-5-fluorophenoxy)-1-piperidinyl]-6-(5-methyl-1,3,4-oxadiazol-2-yl)-pyridazine
MF: C₁₈H₁₇ClFN₅O₂
FW: 389.8
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10566 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10566 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10566 in an organic solvent purged with an inert gas. CAY10566 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10566 in ethanol is approximately 0.15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

CAY10566 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10566 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10566 has a solubility of approximately 0.15 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.

Stearoyl-CoA desaturase (SCD) catalyzes the committed step of the conversion of saturated, long-chain fatty acids to monounsaturated fatty acids. The SCD1 gene is thought to play a key role in lipid homeostasis and body weight regulation.¹⁻³ Thus, modulating SCD1 activity pharmacologically may be a useful tool for regulating type 2 diabetes, dyslipidemia, and obesity. CAY10566 is a potent and selective inhibitor of SCD1 that demonstrates IC₅₀ values of 4.5 and 26 nM in mouse and human enzymatic assays, respectively.⁴ This compound inhibits the conversion of saturated, long-chain fatty acyl-CoAs to monounsaturated, long-chain fatty acyl-CoAs in HepG2 cells with IC₅₀ values of 7.9 and 6.8 nM, respectively, when heptadecanoic acid and palmitic acid are used as the substrate.⁴

References

1. Cohen, P., Miyazaki, M., Socci, N.D., *et al.* Role for stearoyl-CoA desaturase-1 in leptin-mediated weight loss. *Science* **297**, 240-243 (2002).
2. Dobrzyn, P., Sampath, H., Dobrzyn, A., *et al.* Loss of stearoyl-CoA desaturase 1 inhibits fatty acid oxidation and increases glucose utilization in the heart. *Am. J. Physiol. Endocrinol. Metab.* **294**, E357-E364 (2008).
3. Miyazaki, M., Flowers, M.T., Sampath, H., *et al.* Hepatic stearoyl-CoA desaturase-1 deficiency protects mice from carbohydrate-induced adiposity and hepatic steatosis. *Cell Metabolism* **6**, 484-496 (2007).
4. Liu, G., Lynch, J.K., Freeman, J., *et al.* Discovery of potent, selective, orally bioavailable stearoyl-CoA desaturase 1 inhibitors. *J. Med. Chem.* **50**, 3086-3100 (2007).

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

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

MATERIAL 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 Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at 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, 12/05/2011