

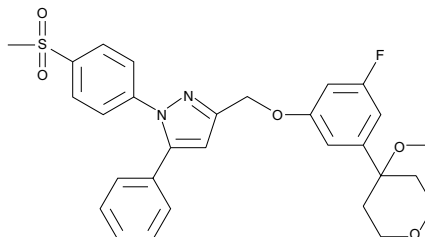
Product Information



CAY10416

Item No. 70635

CAS Registry No.: 443919-96-8
Formal Name: 3-[[[3-fluoro-5-(tetrahydro-4-methoxy-2H-pyran-4-yl)phenoxy]methyl]-1-[4-(methylsulfonyl)phenyl]-5-phenyl-1H-pyrazole
MF: C₂₉H₂₉FN₂O₅S
FW: 536.6
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 204, 269 nm



Laboratory Procedures

For long term storage, we suggest that CAY10416 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10416 is supplied as a crystalline solid. A stock solution may be made by dissolving CAY10416 in an organic solvent purged with an inert gas. CAY10416 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10416 in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Dual cyclooxygenase-2 (COX-2)/5-lipoxygenase (5-LO) inhibitors are potential therapeutic agents for inflammatory diseases and prostate cancer. CAY10416 is a dual COX-2/5-LO inhibitor with IC₅₀ values of 50 and 3 nM, respectively.¹ The selectivity of CAY10416 for COX is greater than 200-fold for COX-2 versus COX-1. COX-2/5-LO inhibitors such as CAY10416 are also apoptosis-inducing agents and are potentially useful in prostate cancer chemotherapy. In the PC3 human carcinoma cell line, CAY10416 exhibits a 45% inhibition of proliferation at 100 μM. In the LNCaP human carcinoma cell line, CAY10416 inhibits growth with an IC₅₀ value of 83 μM.

Reference

1. Barbey, S., Goossens, L., Taverne, T., *et al.* Synthesis and activity of a new methoxytetrahydropyran derivative as dual cyclooxygenase-2/5-lipoxygenase inhibitor. *Bioorg. Medicinal Chem. Letters* **12**, 779-782 (2002).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/70635

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 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, 06/19/2012

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