

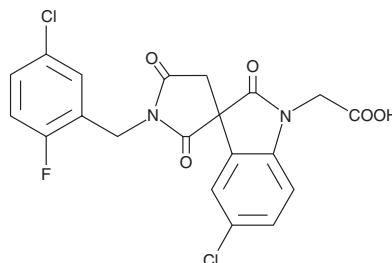
Product Information



CAY10595

Item No. 10012553

CAS Registry No.: 916047-16-0
Formal Name: 5-chloro-1'-[5-chloro-2-fluorophenyl)methyl]-2,2',5'-trioxo-spiro[3H-indole-3,3'pyrrolidine]-1(2H)-acetic acid
MF: C₂₀H₁₃Cl₂FN₂O₅
FW: 451.2
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10595 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10595 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10595 in an organic solvent purged with an inert gas. CAY10595 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10595 in these solvents is approximately 25 mg/ml.

CAY10595 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10595 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10595 has a solubility of approximately 0.5 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

The biological effects of prostaglandin D₂ (PGD₂) are transduced by at least two 7-transmembrane G protein-coupled receptors, designated DP₁ and CRTH2/DP₂. In humans, CRTH2/DP₂ is expressed on Th2 cells, eosinophils, and basophils where it mediates the chemotactic activity of PGD₂.¹ CAY10595, as a racemic mixture, is a potent CRTH2/DP₂ receptor antagonist that binds to the human receptor with a K_i value of 10 nM.² The R-enantiomer of CAY10595 is significantly more potent, exhibiting K_i values of 5.3 and 5 nM at the human and murine CRTH2/DP₂ receptors, respectively. The R-enantiomer of CAY10595 inhibits eosinophil chemotaxis induced by 13,14-dihydro-15-keto-PGD₂ with an IC₅₀ value of 7.3 nM.²

References

1. Hirai, H., Tanaka, K., Yoshie, O., *et al.* Prostaglandin D₂ selectivity induces chemotaxis in T helper type 2 cells, eosinophils, and basophils *via* seven-transmembrane receptor CRTH2. *J. Exp. Med.* **193**(2), 255-261 (2001).
2. Crosignani, S., Page, P., Missotten, M., *et al.* Discovery of a new class of potent, selective, and orally bioavailable CRTH2 (DP₂) receptor antagonists for the treatment of allergic inflammatory diseases. *J. Med. Chem.* **51**, 2227-2243 (2008).

Related Products

15(R)-Prostaglandin D₂ - Item No. 10118 • BAY-u3405 - Item No. 10156 • 11-deoxy-11-methylene-15-keto Prostaglandin D₂ - Item No. 12415 • CAY10648 - Item No. 9000549 • CAY10471 - Item No. 10006735 • CAY10597 - Item No. 10012539

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, 03/22/2011

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