

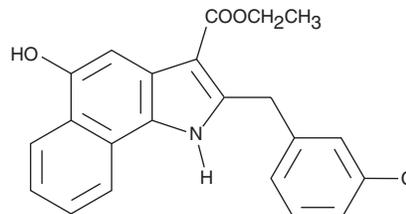
PRODUCT INFORMATION



CAY10606

Item No. 13381

CAS Registry No.: 1159576-98-3
Formal Name: 2-[(3-chlorophenyl)methyl]-5-hydroxy-1H-benz[*g*]indole-3-carboxylic acid, ethyl ester
MF: C₂₂H₁₈ClNO₃
FW: 379.8
Purity: ≥98%
UV/Vis.: λ_{max}: 265, 348 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CAY10606 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10606 in the solvent of choice, which should be purged with an inert gas. CAY10606 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10606 in ethanol is approximately 20 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

5-Lipoxygenase (5-LO) initiates the synthesis of leukotrienes (LTs) from arachidonic acid, primarily in certain leukocyte populations. CAY10606 is a potent, reversible inhibitor of 5-LO, both in cell-free assays (IC₅₀ = 86 nM) and in intact neutrophils (IC₅₀ = 230 nM).¹ It prevents the production of LTs in whole blood, whether 5-LO is activated with the calcium ionophore A23187 (IC₅₀ = 1.6 μM) or formyl peptide (fMLP) following priming with lipopolysaccharide (IC₅₀ = 830 nM). In rats subjected to carrageenan-induced pleurisy, CAY10606 significantly reduces both LT biosynthesis and the inflammatory reaction.¹

Reference

1. Karg, E.M., Luderer, S., Pergola, C., *et al.* Structural optimization and biological evaluation of 2-substituted 5-hydroxyindole-3-carboxylates as potent inhibitors of human 5-lipoxygenase. *J. Med. Chem.* **52**(11), 3474-3483 (2009).

WARNING

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

SAFETY 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, 03/20/2024

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