

PRODUCT INFORMATION

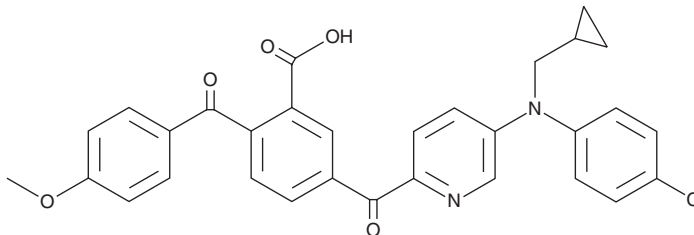


TK05

Item No. 19793

CAS Registry No.: 1245734-61-5
Formal Name: 5-[[5-[(4-chlorophenyl)(cyclopropylmethyl)amino]-2-pyridinyl]carbonyl]-2-(4-methoxybenzoyl)-benzoic acid

MF: $C_{31}H_{25}ClN_2O_5$
FW: 541.0
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Description

TK05 is an inhibitor of leukotriene C_4 (LTC_4) synthase ($IC_{50} = 95$ nM).¹ It is selective for LTC_4 synthase over microsomal prostaglandin E synthase-1 (mPGES-1), microsomal glutathione S-transferase 2 (GST2), and 5-lipoxygenase (5-LO) at 10 μM . TK05 (1 μM) inhibits LTC_4 synthase in Mono-Mac-6 leukemia cells and isolated human eosinophils, monocytes, and platelets but not cord blood-derived mast cells. *In vivo*, TK05 (0.12 and 6 mg/kg) prevents zymosan A-induced increases in peritoneal fluid levels of LTE_4 and LTB_4 in mice.

Reference

1. Kleinschmidt, T.K., Haraldsson, M., Basavarajappa, D., *et al.* Tandem benzophenone amino pyridines, potent and selective inhibitors of human leukotriene C_4 synthase. *J. Pharmacol. Exp. Ther.* **355**(1), 108-116 (2015).

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, 01/30/2024

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