

PRODUCT INFORMATION

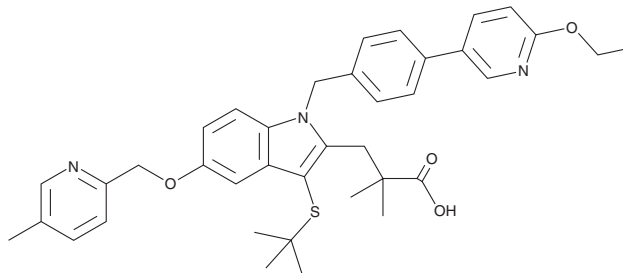


GSK2190915

Item No. 28600

CAS Registry No.: 936350-00-4
Formal Name: 3-[(1,1-dimethylethyl)thio]-1-[[4-(6-ethoxy-3-pyridinyl)phenyl]methyl]- α,α -dimethyl-5-[(5-methyl-2-pyridinyl)methoxy]-1H-indole-2-propanoic acid

Synonym: AM803
MF: C₃₈H₄₃N₃O₄S
FW: 637.8
Purity: \geq 98%
UV/Vis.: λ_{max} : 215, 262 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

GSK2190915 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2190915 in the solvent of choice, which should be purged with an inert gas. GSK2190915 is soluble in organic solvents such as methanol and DMSO.

Description

GSK2190915 is a selective inhibitor of human 5-lipoxygenase-activating protein (FLAP; IC₅₀ = 2.9 nM).¹ It is selective for FLAP over 5-, 12-, and 15-lipoxygenase, leukotriene A₄ (LTA₄) hydrolase, LTC₄ synthase, COX-1, and COX-2 (IC₅₀s = >10 μ m for all). GSK2190915 inhibits production of LTB₄ induced by the calcium ionophore A23187 (Item No. 11016) in isolated whole blood from human, rat, and mouse (IC₅₀s = 436, 223, and 148 nM, respectively). It decreases increases in the levels of cysteinyl leukotriene (CysLT) and LTB₄ induced by lung instillation of A23187 in the bronchoalveolar lavage fluid (BALF) of rats when administered at doses of 1 and 3 mg/kg. GSK2190915 (3 mg/kg) also reduces increases in the levels of CysLT, LTB₄, myeloperoxidase (MPO), and plasma protein extravasation in a mouse model of peritonitis induced by zymosan in mice.

Reference

1. Lorrain, D.S., Bain, G., Correa, L.D., *et al.* Pharmacology of AM803, a novel selective five-lipoxygenase-activating protein (FLAP) inhibitor in rodent models of acute inflammation. *Eur. J. Pharmacol.* **640**(1-3), 211-218 (2010).

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, 11/08/2022

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