

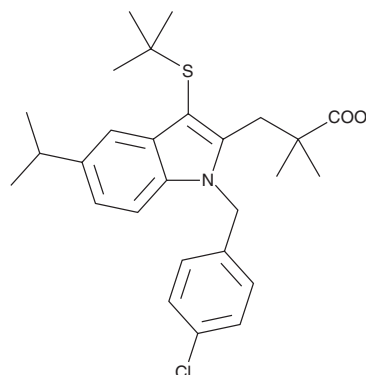
PRODUCT INFORMATION



MK-886

Item No. 21753

CAS Registry No.: 118414-82-7
Formal Name: 1-[(4-chlorophenyl)methyl]-3-[(1,1-dimethylethyl)thio]- α,α -dimethyl-5-(1-methylethyl)-1H-indole-2-propanoic acid
Synonym: L-663,536
MF: C₂₇H₃₄ClNO₂S
FW: 472.1
Purity: \geq 98%
UV/Vis.: λ_{max} : 225, 288 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

MK-886 is supplied as a crystalline solid. A stock solution may be made by dissolving the MK-886 in the solvent of choice. MK-886 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of MK-886 in these solvents is approximately 0.14, 5, and 15 mg/ml, respectively.

MK-886 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MK-886 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MK-886 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MK-886 is a potent, selective inhibitor of 5-lipoxygenase-activating protein (FLAP) that prevents 5-lipoxygenase activation *in vivo* and inhibits leukotriene biosynthesis in leukocytes with an IC₅₀ of 2.5 nM.¹ In human whole blood, leukotriene biosynthesis is inhibited by MK-886 with an IC₅₀ of 1.1 μ M.

Reference

1. Dixon, R.A.F., Diehl, R.E., Opas, E., *et al.* Requirement of a 5-lipoxygenase-activating protein for leukotriene synthesis. *Nature* **343**, 282-284 (1990).

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, 10/31/2022

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