

PRODUCT INFORMATION

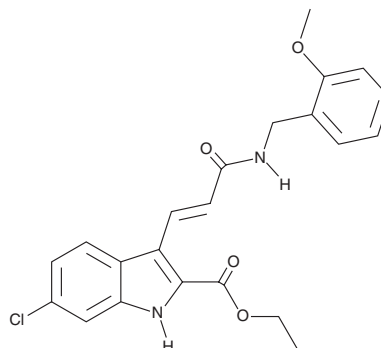


9c(i472)

Item No. 28225

CAS Registry No.: 2349374-37-2
Formal Name: 6-chloro-3-[(1E)-3-[[[2-methoxyphenyl)methyl]amino]-3-oxo-1-propen-1-yl]-1H-indole-2-carboxylic acid, ethyl ester

Synonym: 15-LOX-1 Inhibitor i472
MF: C₂₂H₂₁ClN₂O₄
FW: 412.9
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 240, 283, 328 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

9c(i472) is supplied as a crystalline solid. A stock solution may be made by dissolving the 9c(i472) in the solvent of choice, which should be purged with an inert gas. 9c(i472) is soluble in the organic solvent DMSO. 9c(i472) is also slightly soluble in methanol.

Description

9c(i472) is an inhibitor of 15-lipoxygenase-1 (15-LO-1; IC₅₀ = 0.19 μM).¹ It decreases LPS- and IFN-γ-induced NF-κB activity in RAW-Blue™ cells when used at concentrations of 0.2, 1, and 5 μM. 9c(i472) reduces LPS- and IFN-γ-induced increases in Nos2 expression and lipid peroxidation in RAW 264.7 cells when used at a concentration of 5 μM.

Reference

1. Guo, H., Verhoek, I.C., Prins, G.G.H., et al. Novel 15-lipoxygenase-1 inhibitor protects macrophages from lipopolysaccharide-induced cytotoxicity. *J. Med. Chem.* **62(9)**, 4624-4637 (2019).

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, 03/18/2024

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