

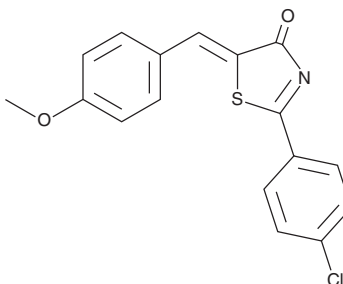
PRODUCT INFORMATION



CAY10649

Item No. 10804

CAS Registry No.: 1272519-89-7
Formal Name: (Z)-2-(4-chlorophenyl)-5-(4-methoxybenzylidene)thiazol-4(5H)-one
MF: C₁₇H₁₂ClNO₂S
FW: 329.8
Purity: ≥98%
UV/Vis.: λ_{max}: 297, 428 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

CAY10649 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10649 in the solvent of choice, which should be purged with an inert gas. CAY10649 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10649 in these solvents is approximately 0.5 mg/ml and 1 mg/ml, respectively.

Description

5-Lipoxygenase (5-LO) catalyzes the biosynthesis of leukotrienes, which are involved in a variety of inflammatory responses, including neutrophil chemotaxis, vascular permeability, and smooth muscle contraction.¹ CAY10649 is a thiazolinone compound that demonstrates direct inhibition of 5-LO product formation in intact polymorphonuclear leukocytes (PMNL) (IC₅₀ = 0.28 μM) and a soluble fraction of a S100 PMNL cell lysate (IC₅₀ = 0.09 μM) after stimulation with calcium and arachidonic acid.²

References

1. Samuelsson, B., Dahlén, S.-E., Lindgren, J.Å., *et al.* Leukotrienes and lipoxins: Structures, biosynthesis, and biological effects. *Science* **237**(4819), 1171-1176 (1987).
2. Hoffman, B., Barzen, S., Rödl, C.B., *et al.* A class of 5-benzylidene-2-phenylthiazolinones with high potency as direct 5-lipoxygenase inhibitors. *J. Med. Chem.* **54**(6), 1943-1947 (2011).

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, 02/21/2024

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