

PRODUCT INFORMATION



ZLJ-6

Item No. 13271

CAS Registry No.: 1051931-39-5
Formal Name: (5Z)-2-amino-1,5-dihydro-1-methyl-5-[[4-(methylsulfonyl)phenyl]methylene]-4H-imidazol-4-one, monomethanesulfonate

MF: C₁₂H₁₃N₃O₃S • CH₃SO₃H
FW: 375.4

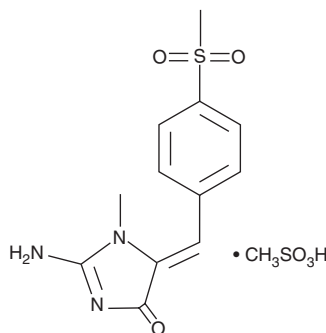
Purity: ≥98%

UV/Vis.: λ_{max}: 240, 338 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

ZLJ-6 is supplied as a crystalline solid. A stock solution may be made by dissolving the ZLJ-6 in an organic solvent purged with an inert gas. ZLJ-6 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ZLJ-6 in these solvents is approximately 10 mg/ml.

ZLJ-6 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ZLJ-6 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ZLJ-6 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ZLJ-6 is a dual inhibitor of COX and 5-lipoxygenase (5-LO) (IC₅₀ = 0.73, 0.31, and 0.99 μM for COX-1, COX-2, and 5-LO, respectively, in whole blood).¹ It reduces carrageenin-induced paw edema (inflammation) in rats and acetic acid-induced abdominal constriction (pain) in mice without causing gastrointestinal ulcers, a common side effect of non-selective COX inhibitors.

Reference

1. Li, L., Ji, H., Sheng, L., *et al.* The anti-inflammatory effects of ZLJ-6, a novel dual cyclooxygenase/5-lipoxygenase inhibitor. *Eur. J. Pharmacol.* **607**(1-3), 244-250 (2009).

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, 03/14/2024

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