

PRODUCT INFORMATION

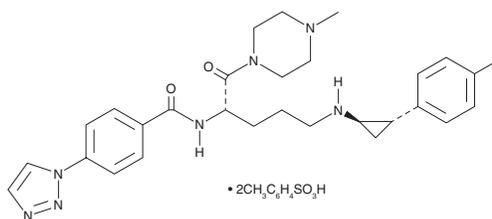


IMG-7289 (tosylate)

Item No. 43733

CAS Registry No.: 1990504-72-7
Formal Name: N-[(1S)-4-[[[(1R,2S)-2-(4-fluorophenyl)cyclopropyl]amino]-1-[4-methyl-1-piperazinyl]carbonyl]butyl]-4-(1H-1,2,3-triazol-1-yl)-benzamide, 4-methylbenzenesulfonate (1:2)

Synonym: Bomedemstat ditosylate
MF: C₂₈H₃₄FN₇O₂ • 2C₇H₈O₃S
FW: 864.0
Purity: ≥98%
Supplied as: A 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

IMG-7289 (tosylate) is supplied as a solid. A stock solution may be made by dissolving the IMG-7289 (tosylate) in the solvent of choice, which should be purged with an inert gas. IMG-7289 (tosylate) is soluble (≥10 mg/ml) in DMSO.

Description

IMG-7289 is an irreversible inhibitor of lysine-specific demethylase 1 (LSD1).¹ It induces cell cycle arrest at the S phase in SET-2 acute megakaryoblastic leukemia cells and apoptosis in Ba/F3 murine hematopoietic cells expressing Jak2 containing a valine-to-phenylalanine substitution at position 617 (Jak2^{V617F}) when used at a concentration of 1 μM. IMG-7289 (45 mg/kg per day) decreases blood levels of leukocytes, neutrophils, monocytes, and platelets, increases bone marrow cell levels of trimethylated lysine 4 histone 3 (H3K4me3) and H3K9me2, decreases spleen fibrosis and weight, and increases survival in a Jak2^{V617F} mouse model of myelofibrosis.

Reference

1. Jutzi, J.S., Kleppe, M., Dias, J., *et al.* LSD1 inhibition prolongs survival in mouse models of MPN by selectively targeting the disease clone. *Hemisphere* **2**(3), e54 (2018).

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, 08/08/2025

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