

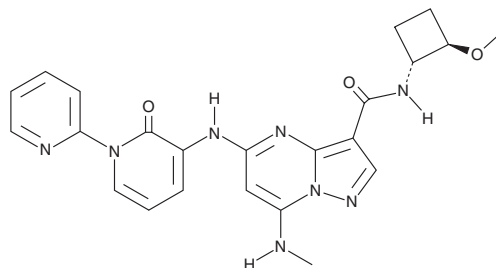
PRODUCT INFORMATION



Zasocitinib

Item No. 35682

CAS Registry No.: 2272904-53-5
Formal Name: N-[(1R,2R)-2-methoxycyclobutyl]-7-(methylamino)-5-[(2-oxo[1(2H),2'-bipyridin]-3-yl)amino]-pyrazolo[1,5-a]pyrimidine-3-carboxamide
Synonyms: NDI-034858, TAK-279
MF: C₂₃H₂₄N₈O₃
FW: 460.5
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

Zasocitinib is supplied as a solid. A stock solution may be made by dissolving the zasocitinib in the solvent of choice, which should be purged with an inert gas. Zasocitinib is soluble in acetonitrile and DMSO.

Description

Zasocitinib is an inhibitor of tyrosine kinase 2 (TYK2).¹ It selectively binds to the TYK2 Janus homology (JH2) domain ($K_d = 0.0038$ nM) over the JAK1 and JAK2 JH2 domains (K_d s = 4,975 and 23,000 nM, respectively). Zasocitinib inhibits the production of chemokine (C-X-C motif) ligand 10 (CXCL10) induced by IFN- α in isolated human, mouse, and rat whole blood (IC_{50} s = 22, 347, and 91 nM, respectively). It also inhibits IL-23-induced STAT3 phosphorylation in human CD161⁺CD3⁺ Th17 cells ($IC_{50} = 3.7$ nM). Zasocitinib (3, 10, and 30 mg/kg twice per day) decreases ankle diameter in a rat model of adjuvant-induced arthritis. It reduces the colon weight-to-length ratio and improves histology scores in a mouse model of colitis induced by an anti-CD40 antibody when administered at doses of 30 and 90 mg/kg.

Reference

1. Leit, S., Greenwood, J., Carriero, S., *et al.* Discovery of a potent and selective tyrosine kinase 2 inhibitor: TAK-279. *J. Med. Chem.* **66**(15), 10473-10496 (2023).

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, 08/20/2024

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