

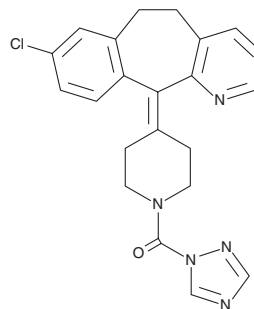
PRODUCT INFORMATION



JZP 361

Item No. 43994

CAS Registry No.: 1680193-80-9
Formal Name: [4-(8-chloro-5,6-dihydro-11H-benzo[5,6]cyclohepta[1,2-b]pyridin-11-ylidene)-1-piperidinyl]-1H-1,2,4-triazol-1-yl-methanone
MF: C₂₂H₂₀ClN₅O
FW: 405.9
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

JZP 361 is supplied as a solid. A stock solution may be made by dissolving the JZP 361 in the solvent of choice, which should be purged with an inert gas. JZP 361 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

JZP 361 is an inhibitor of monoacylglycerol lipase (MAGL; IC₅₀ = 46 nM).¹ It is selective for MAGL over fatty acid amide hydrolase (FAAH) and α/β-hydrolase domain-containing protein 6 (ABHD6; IC₅₀s = 7,240 and 1,790 nM, respectively). JZP 361 also antagonizes histamine activity in isolated guinea pig ileum (pA₂ = 6.81).

Reference

1. Patel, J.Z., Ahenkorah, S., Vaara, M., *et al.* Loratadine analogues as MAGL inhibitors. *Bioorg. Med. Chem. Lett.* **25**(7), 1436-1442 (2015).

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, 08/26/2025

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