

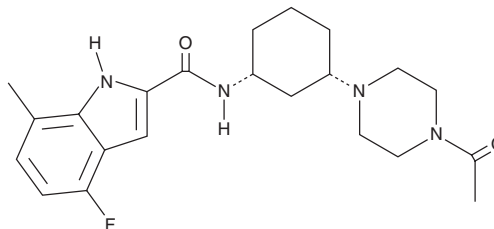
PRODUCT INFORMATION



EZM0414

Item No. 43914

CAS Registry No.: 2411748-50-8
Formal Name: N-[(1R,3S)-3-(4-acetyl-1-piperazinyl)cyclohexyl]-4-fluoro-7-methyl-1H-indole-2-carboxamide
MF: C₂₂H₂₉FN₄O₂
FW: 400.5
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

EZM0414 is supplied as a solid. A stock solution may be made by dissolving the EZM0414 in the solvent of choice, which should be purged with an inert gas. EZM0414 is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

EZM0414 is an inhibitor of SET domain-containing protein 2 (SETD2; IC₅₀ = 18 nM).¹ It inhibits the proliferation of A549 non-small cell lung cancer (NSCLC) cells (EC₅₀ = 57 nM). EZM0414 (15 and 30 mg/kg) decreases tumor volume in a KMS-11 multiple myeloma mouse xenograft model. It decreases cardiomyocyte levels of trimethylated lysine 36 on histone H3 (H3K36Me3) in a mouse model of heart failure with preserved ejection fraction (HFpEF) induced by a high-fat diet and L-NAME (Item No. 80210).² EZM0414 also reduces lung congestion and cardiac hypertrophy in the same model.

References

1. Alford, J.S., Lampe, J.W., Brach, D., *et al.* Conformational-design-driven discovery of EZM0414: A selective, potent SETD2 inhibitor for clinical studies. *ACS Med. Chem. Lett.* **13**(7), 1137-1143 (2022).
2. Costantino, S., Mohammed, S.A., Ambrosini, S., *et al.* Chromatin rewiring by SETD2 drives lipotoxic injury in cardiometabolic HFpEF. *Circ. Res.* **136**(10), 1079-1095 (2025).

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, 09/17/2025

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