

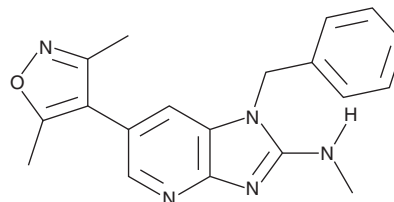
PRODUCT INFORMATION



ZEN-3694

Item No. 43661

CAS Registry No.: 1643947-30-1
Formal Name: 6-(3,5-dimethyl-4-isoxazolyl)-N-methyl-1-(phenylmethyl)-1H-imidazo[4,5-b]pyridin-2-amine
MF: C₁₉H₁₉N₅O
FW: 333.4
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

ZEN-3694 is supplied as a solid. A stock solution may be made by dissolving the ZEN-3694 in the solvent of choice, which should be purged with an inert gas. ZEN-3694 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

ZEN-3694 is an inhibitor of the bromodomain and extra terminal domain (BET) family of proteins.^{1,2} It induces apoptosis in wild-type MCF-7 cells and those resistant to palbociclib (PD 0332991; Item No. 16273) when used at a concentration of 1 μM.²

References

1. Feng, Y., Mahdi, H., Piekarz, R., *et al.* An LC-MS/MS method for determination of the bromodomain inhibitor ZEN-3694 and its metabolite ZEN-3791 in human plasma. *Bioanalysis* **16**(8), 227-238 (2024).
2. Kharenko, O.A., Patel, R.G., Calosing, C., *et al.* Combination of ZEN-3694 with CDK4/6 inhibitors reverses acquired resistance to CDK4/6 inhibitors in ER-positive breast cancer. *Cancer Gene Ther.* **29**(6), 859-869 (2022).

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, 06/03/2025

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