

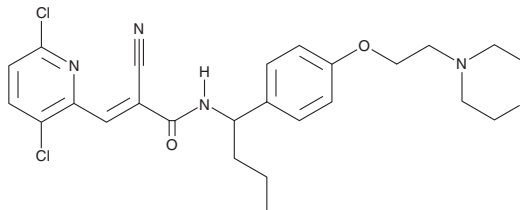
PRODUCT INFORMATION



EOAI3402143

Item No. 43117

CAS Registry No.: 1699750-95-2
Formal Name: 2E-cyano-3-(3,6-dichloro-2-pyridinyl)-
N-[1-[4-[2-(4-morpholinyl)ethoxy]
phenyl]butyl]-2-propenamide
Synonym: G9
MF: C₂₅H₂₈Cl₂N₄O₃
FW: 503.4
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

EOAI3402143 is supplied as a solid. A stock solution may be made by dissolving the EOAI3402143 in the solvent of choice, which should be purged with an inert gas. EOAI3402143 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

EOAI3402143 is an inhibitor of ubiquitin-specific protease 9X (USP9X), USP24, and USP5.¹ It inhibits the catalytic domain of USP9X (IC₅₀ = 1.6 μM). EOAI3402143 (5 μM) inhibits USP9X, USP24, and USP5 in MM.1S multiple myeloma cells. It inhibits the growth of vemurafenib-sensitive and -resistant A375 melanoma cells (IC₅₀s = 1 μM).² *In vivo*, EOAI3402143 (15 mg/kg) reduces tumor volume in MM.1S and A375 mouse xenograft models.^{1,2}

References

1. Peterson, L.F., Sun, H., Liu, Y., *et al.* Targeting deubiquitinase activity with a novel small-molecule inhibitor as therapy for B-cell malignancies. *Blood* **125**(23), 3588-3597 (2015).
2. Potu, H., Peterson, L.F., Pal, A., *et al.* Usp5 links suppression of p53 and FAS levels in melanoma to the BRAF pathway. *Oncotarget* **5**(14), 5559-5569 (2014).

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, 04/08/2025

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