

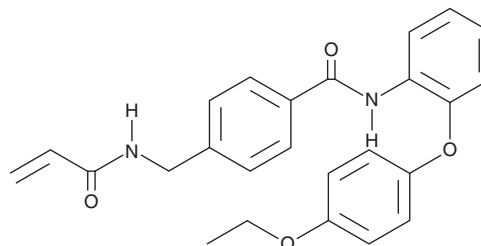
PRODUCT INFORMATION



EN4

Item No. 33473

CAS Registry No.: 1197824-15-9
Formal Name: N-[2-(4-ethoxyphenoxy)phenyl]-4-[[[(1-oxo-2-propen-1-yl)amino]methyl]-benzamide
MF: C₂₅H₂₄N₂O₄
FW: 416.5
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

EN4 is supplied as a solid. A stock solution may be made by dissolving the EN4 in the solvent of choice, which should be purged with an inert gas. EN4 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

EN4 is a covalent inhibitor of c-Myc.¹ It inhibits c-Myc-induced increases in transcriptional activity in a reporter assay using HEK293T cells (EC₅₀ = 12.5 μM). EN4 (1, 10, or 50 μM) reduces the proliferation of 231MFP breast cancer cells. *In vivo*, EN4 (50 mg/kg per day) decreases tumor volume in a 231MFP mouse xenograft model.

Reference

1. Boike, L., Cioffi, A.G., Majewski, F.C., *et al.* Discovery of a functional covalent ligand targeting an intrinsically disordered cysteine within MYC. *Cell Chem. Biol.* **28(1)**, 4-13 (2021).

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, 10/22/2024

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