PRODUCT INFORMATION



AOH1996

Item No. 39526

CAS Registry No.: Formal Name:	2089314-64-5 N-[2-[[2-(3-methoxyphenoxy)phenyl]amino]-	
	2-oxoethyij-1-haphthalehecarboxamide	
MF:	$C_{26}H_{22}N_{2}O_{4}$	
FW:	426.5	N H
Purity:	≥98%	
UV/Vis.:	λ _{max} : 222 nm	
Supplied as:	A solid	
Storage:	-20°C	$\langle \rangle \rangle$
Stability:	≥4 years	
Information represents the product specifications. Patch specific analytical results are provided on each certificate of analysis		

Laboratory Procedures

AOH1996 is supplied as a solid. A stock solution may be made by dissolving the AOH1996 in the solvent of choice, which should be purged with an inert gas. AOH1996 is soluble in ethanol, methanol, and chloroform.

Description

AOH1996 is an inhibitor of proliferating cell nuclear antigen (PCNA).¹ It binds to PCNA in a thermal shift assay. AOH1996 (500 nM) induces apoptosis in SK-N-DZ neuroblastoma cells but not 7SMOO32 normal neural crest stem cells. It induces DNA damage in SK-N-AS neuroblastoma cells expressing wild-type DNA-directed RNA polymerase II subunit RPB1 but not in SK-N-AS cells expressing RPB1^{Y418A}. AOH1996 (40 mg/kg) decreases tumor volume in SK-N-BE(2)-C neuroblastoma, MDA-MB-468 breast cancer, and NCI H82 small cell lung cancer mouse xenograft models.

Reference

1. Gu, L., Li, M., Li, C.M., et al. Small molecule targeting of transcription-replication conflict for selective chemotherapy. Cell Chem. Biol. 30, 1-13 (2023).

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.

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, 02/24/2025

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