

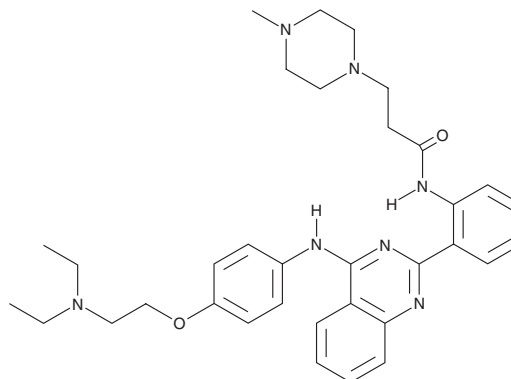
PRODUCT INFORMATION



hVEGF-IN-1

Item No. 37550

CAS Registry No.: 1637443-98-1
Formal Name: N-[2-[4-[[4-[2-(diethylamino)ethoxy]phenyl]amino]-2-quinazoliny]phenyl]-4-methyl-1-piperazinepropanamide
Synonym: Human Vascular Endothelial Growth Factor Inhibitor 1
MF: C₃₄H₄₃N₇O₂
FW: 581.8
Purity: ≥98%
UV/Vis.: λ_{max}: 241 nm
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

hVEGF-IN-1 is supplied as a solid. A stock solution may be made by dissolving the hVEGF-IN-1 in the solvent of choice, which should be purged with an inert gas. hVEGF-IN-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of hVEGF-IN-1 in DMSO is approximately 2 mg/ml and approximately 25 mg/ml in ethanol and DMF.

Description

hVEGF-IN-1 is an inhibitor of VEGFA translation.¹ It binds to the mRNA encoding VEGF-A ($K_d = 4.64 \mu\text{M}$) and decreases its translation when used at a concentration of 1.5 μM . hVEGF-IN-1 (3 μM) reduces migration of MCF-7 breast cancer cells. It decreases tumor growth and weight without reducing body weight in an MCF-7 mouse xenograft model when administered at a dose of 7.5 mg/kg per day.

Reference

1. Wang, S.-K., Wu, Y., Wang, X.-Q., *et al.* Discovery of small molecules for repressing cap-independent translation of human vascular endothelial growth factor (hVEGF) as novel antitumor agents. *J. Med. Chem.* **60**(13), 5306-5319 (2017).

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, 06/09/2023

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