

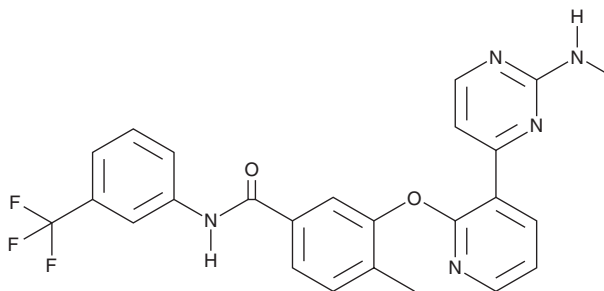
PRODUCT INFORMATION



AMG-Tie2-1

Item No. 28850

CAS Registry No.: 870223-96-4
Formal Name: 4-methyl-3-[[3-[2-(methylamino)-4-pyrimidinyl]-2-pyridinyl]oxy]-N-[3-(trifluoromethyl)phenyl]-benzamide
MF: C₂₅H₂₀F₃N₅O₂
FW: 479.5
Purity: ≥98%
UV/Vis.: λ_{max}: 245 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

AMG-Tie2-1 is supplied as a solid. A stock solution may be made by dissolving the AMG-Tie2-1 in the solvent of choice, which should be purged with an inert gas. AMG-Tie2-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AMG-Tie2-1 in these solvents is approximately 10 mg/ml.

AMG-Tie2-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AMG-Tie2-1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AMG-Tie2-1 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AMG-Tie2-1 is an inhibitor of tunica interna endothelial cell kinase 2 (Tie2) and VEGF receptor 2 (VEGFR2) with IC₅₀ values of 1 and 3 nM, respectively, in a homogenous time-resolved fluorescence (HTRF) assay.¹ It inhibits angiotensin 1-induced Tie2 autophosphorylation in EA.hy926 human endothelial cells (IC₅₀ = 10 nM).

Reference

1. Hodous, B.L., Geuns-Meyer, S.D., Hughes, P.E., *et al.* Evolution of a highly selective and potent 2-(pyridin-2-yl)-1,3,5-triazine Tie-2 kinase inhibitor. *J. Med. Chem.* **50**(4), 611-626 (2007).

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, 11/23/2022

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