

PRODUCT INFORMATION

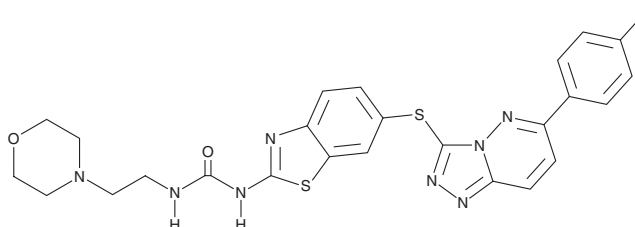


SAR125844

Item No. 37542

CAS Registry No.: 1116743-46-4
Formal Name: N-[6-[[6-(4-fluorophenyl)-1,2,4-triazolo[4,3-b]pyridazin-3-yl]thio]-2-benzothiazolyl]-N'-[2-(4-morpholinyl)ethyl]-urea

MF: C₂₅H₂₃FN₈O₂S₂
FW: 550.6
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 282 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

SAR125844 is supplied as a solid. A stock solution may be made by dissolving the SAR125844 in the solvent of choice, which should be purged with an inert gas. SAR125844 is slightly soluble in DMSO and dimethyl formamide.

Description

SAR125844 is an inhibitor of the receptor tyrosine kinase c-Met (IC₅₀ = 4 nM).¹ It is selective for c-Met over tropomyosin-related kinase A (TrkA), TrkB, PDGFRα, Axl, and Mer (IC₅₀s = 39, 280, 55, 87, and 105 nM, respectively).² SAR125844 inhibits the proliferation of MKN45 gastric cancer cells (IC₅₀ = 7 nM).¹ *In vivo*, SAR125844 (20 mg/kg) inhibits intratumoral c-Met phosphorylation in a MET-amplified Hs 746T gastric cancer mouse xenograft model. SAR125844 (10-45 mg/kg) also reduces tumor volume in Hs 746T and SNU-5 gastric cancer mouse xenograft models.²

References

1. Ugolini, A., Kenigsberg, M., Rak, A., *et al.* Discovery and pharmacokinetic and pharmacological properties of the potent and selective MET kinase inhibitor 1-{6-[[6-(4-Fluorophenyl)-[1,2,4]triazolo[4,3-b]pyridazin-3-ylsulfanyl]benzothiazol-2-yl]-3-(2-morpholin-4-ylethyl)urea (SAR125844). *J. Med. Chem.* **59(15)**, 7066-7074 (2016).
2. Egile, C., Kenigsberg, M., Delaisi, C., *et al.* The selective intravenous inhibitor of the MET tyrosine kinase SAR125844 inhibits tumor growth in MET-amplified cancer. *Mol. Cancer Ther.* **14(2)**, 384-394 (2015).

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, 04/28/2023

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