

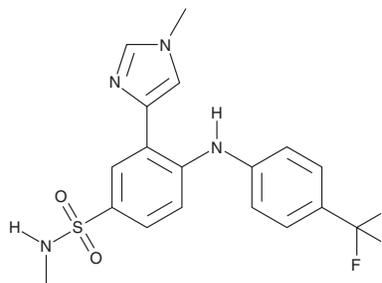
PRODUCT INFORMATION



VT103

Item No. 42327

CAS Registry No.: 2290608-13-6
Formal Name: N-methyl-3-(1-methyl-1H-imidazol-4-yl)-4-[[4-(trifluoromethyl)phenyl]amino]benzenesulfonamide
MF: C₁₈H₁₇F₃N₄O₂S
FW: 410.4
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

VT103 is supplied as a solid. A stock solution may be made by dissolving the VT103 in the solvent of choice, which should be purged with an inert gas. VT103 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

VT103 is an inhibitor of TEA domain transcription factor 1 (TEAD1; IC₅₀ = 1.02 nM).¹ It inhibits TEAD autopalmitoylation in HEK293 cells expressing TEAD1 but not HEK293 cells expressing TEAD2, TEAD3, or TEAD4 when used at a concentration of 3 μM. VT103 also inhibits palmitoylation of endogenous TEAD1 in NCI-H2373 pleural mesothelioma cells. *In vivo*, VT103 (0.3-10 mg/kg) reduces tumor volume in NCI-H2373 and NCI H226 mesothelioma mouse xenograft models.

Reference

1. Tang, T.T., Konradi, A.W., Feng, Y., *et al.* Small molecule inhibitors of TEAD auto-palmitoylation selectively inhibit proliferation and tumor growth of NF2-deficient mesothelioma. *Mol. Cancer Ther.* **20(6)**, 986-998 (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, 12/17/2024

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