

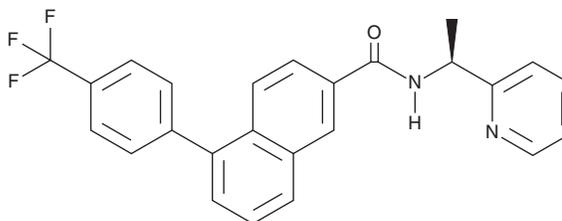
# PRODUCT INFORMATION



VT104

Item No. 45223

CAS Registry No.: 2417718-25-1  
Formal Name: N-[(1S)-1-(2-pyridinyl)ethyl]-5-[4-(trifluoromethyl)phenyl]-2-naphthalenecarboxamide  
MF: C<sub>25</sub>H<sub>19</sub>F<sub>3</sub>N<sub>2</sub>O  
FW: 420.4  
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

VT104 is supplied as a solid. A stock solution may be made by dissolving the VT104 in the solvent of choice, which should be purged with an inert gas. VT104 is soluble (≥10 mg/ml) in ethanol and DMSO.

## Description

VT104 is an inhibitor of TEA domain transcription factor (TEAD) autopalmitoylation.<sup>1</sup> It inhibits autopalmitoylation of TEAD1-4 in HEK293 cells expressing the human transcription factors when used at a concentration of 3 μM. VT104 inhibits the growth of NCI-H2052, NCI-H2373, and NCI H226 mesothelioma cells (IC<sub>50</sub>s = 31.6, 25.6, and 16.1 nM, respectively). *In vivo*, VT104 (1, 3, and 10 mg/kg) reduces tumor volume in an NCI H226 mouse xenograft model.

## 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, 02/23/2026

## 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