

PRODUCT INFORMATION



YAP/TAZ Inhibitor-3

Item No. 45146

CAS Registry No.: 2506273-81-8
Formal Name: N-[(1R)-2-hydroxy-1-methylethyl]-5-[4-(trifluoromethyl)phenoxy]-2-naphthalenecarboxamide

Synonyms: VT3989, Yes-associated Protein/Transcriptional Coactivator with PDZ-binding Domain Inhibitor-3

MF: C₂₁H₁₈F₃NO₃

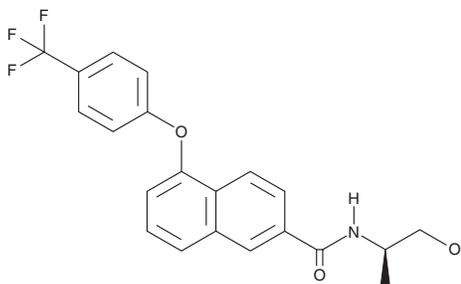
FW: 389.4

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

YAP/TAZ inhibitor-3 is supplied as a solid. A stock solution may be made by dissolving the YAP/TAZ inhibitor-3 in the solvent of choice, which should be purged with an inert gas. YAP/TAZ inhibitor-3 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

YAP/TAZ inhibitor-3 is an inhibitor of the protein-protein interaction between TEA domain transcription factor (TEAD) and the transcriptional coactivators yes-associated transcriptional regulator (YAP) and its paralog, transcriptional coactivator with PDZ-binding motif (TAZ).^{1,2} It inhibits TEAD1-4 palmitoylation, which is essential for the TEAD-YAP/TAZ interaction, in HEK293T cells expressing TEAD1-4 when used at a concentration of 3 μM. YAP/TAZ inhibitor-3 inhibits the proliferation of a variety of mesothelioma cells, including those with homozygous *NF2* mutations (GI₅₀s = 8-74 nM) or heterozygous *NF2* truncations (GI₅₀s = 178-1,511 nM), both of which lead to constitutive TEAD activation, as well as some mesothelioma cells expressing wild-type *NF2* (GI₅₀s = 38->3,000 nM).¹ It also reduces tumor growth in an NCI H226 mouse xenograft model when administered at doses of 3 and 10 mg/kg.

References

1. Yap, T.A., Kwiatkowski, D.J., Dagogo-Jack, I., *et al.* YAP/TEAD inhibitor VT3989 in solid tumors: A phase 1/2 trial. *Nat. Med.* **31(12)**, 4281-4290 (2025).
2. Chen, F., Su, J., Liu, Y., *et al.* Targeting YAP/TAZ-TEAD and their protein-protein interaction for precision cancer therapy. *Eur. J. Med. Chem.* **302(Pt. 2)**, (2026).

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, 02/19/2026

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