

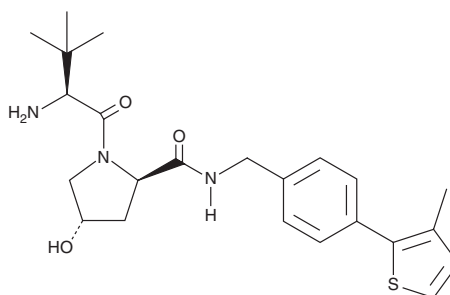
PRODUCT INFORMATION



(R,S,S)-VH 032

Item No. 44838

CAS Registry No.: 2230826-33-0
Formal Name: (4S)-3-methyl-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-D-prolinamide
MF: C₂₂H₃₀N₄O₃S
FW: 430.6
Purity: ≥95%
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

(R,S,S)-VH 032 is supplied as a solid. A stock solution may be made by dissolving the (R,S,S)-VH 032 in the solvent of choice, which should be purged with an inert gas. (R,S,S)-VH 032 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

(R,S,S)-VH 032 is a diastereomer of VHL ligand 1 (Item No. 21591), a building block in the synthesis of proteolysis-targeting chimeras (PROTACs).^{1,2} (R,S,S)-VH 032-containing molecules bind targets with the same affinity as PROTACs utilizing VHL ligand 1 but do not induce target degradation and have been used as controls for VHL ligand 1-containing PROTAC degradation activity.

References

1. Burslem, G.M., Smith, B.E., Lai, A.C., *et al.* The advantages of targeted protein degradation over inhibition: An RTK case study. *Cell Chem. Biol.* **25(1)**, 66-77 (2018).
2. Burslem, G.M., Song, J., Chen, X., *et al.* Enhancing antiproliferative activity and selectivity of a FLT-3 inhibitor by proteolysis targeting chimera conversion. *J. Am. Chem. Soc.* **140(48)**, 16428-16432 (2018).

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, 03/03/2026

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