

PRODUCT INFORMATION

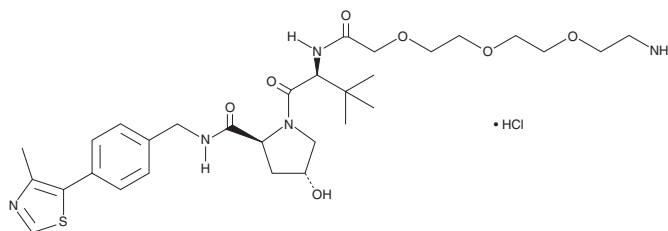


VH 032 Linker 1

Item No. 26148

CAS Registry No.: 2097971-11-2
Formal Name: (4R)-N-[2-[2-[2-(2-aminoethoxy)ethoxy]ethoxy]acetyl]-3-methyl-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide, monohydrochloride

MF: C₃₀H₄₅N₅O₇S • HCl
FW: 656.2
Purity: ≥95%
UV/Vis.: λ_{max}: 272 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

VH 032 linker 1 is a precursor in the synthesis of the proteolysis-targeting chimeras (PROTACs) that contain a terminal amine conjugated via a PEG linker to VHL ligand 1 (Item No. 21591), which binds to VHL E3 ligase. It has been used in the synthesis of PROTACs targeting TEA domain transcription factors (TEADs) in cancer cells.¹

Reference

1. Zbeig, J.R., Villemure, E., Rudolph, J., *et al.* Therapeutic compounds and methods of use. *Genentech, Inc.* WO2021178339AI (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, 01/25/2024

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