

PRODUCT INFORMATION



VH 032 Linker 2

Item No. 26149

CAS Registry No.: 2064292-52-8
Formal Name: (2S,4R)-1-((S)-17-amino-2-(tert-butyl)-4-oxo-6,9,12,15-tetraoxa-3-azaheptadecanoyl)-4-hydroxy-N-(4-(4-methylthiazol-5-yl)benzyl)pyrrolidine-2-carboxamide, monohydrochloride

Synonym: E3 Ligase Ligand-Linker Conjugate 7

MF: C₃₂H₄₉N₅O₈S • HCl

FW: 700.3

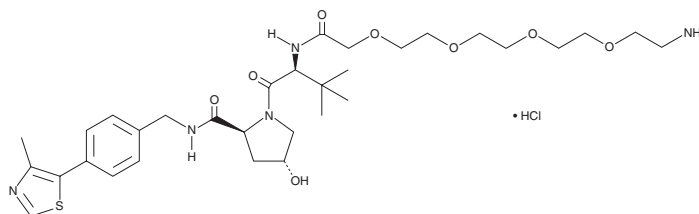
Purity: ≥98%

UV/Vis.: λ_{max}: 271 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

VH 032 Linker 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the VH 032 linker 2 in the solvent of choice. VH 032 linker 2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of VH 032 linker 2 in ethanol is approximately 0.7 mg/ml and 1.1 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of VH 032 linker 2 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of VH 032 linker 2 in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

VH 032 Linker 2 is a derivative of the proteolysis-targeting chimera technology (PROTAC) building block VHL ligand 1 (Item No. 21591). VH 032 Linker 2 comprises the von Hippel-Lindau (VHL) ligand domain and the polyethylene glycol (PEG) linker of the bromodomain and extra terminal (BET) inhibitor-containing PROTACs MZ2 and MZP-55.¹

Reference

1. Chan, K.H., Zengerle, M., and Ciulli, A. Impact of target warhead and linkage vector on inducing protein degradation: Comparison of bromodomain and extra-terminal (BET) degraders derived from triazolodiazepine (JQ1) and tetrahydroquinoline (l-BET726) BET inhibitor scaffolds. *J. Med. Chem.* **61**(2), 504-513 (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, 01/21/2019

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