

PRODUCT INFORMATION



VH 032 Linker 3 (hydrate)

Item No. 26151

Formal Name: (4R)-N-(5-amino-1-oxopentyl)-3-methyl-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide, dihydrochloride, hydrate

MF: C₂₇H₃₉N₅O₄S • 2HCl [XH₂O]

FW: 602.6

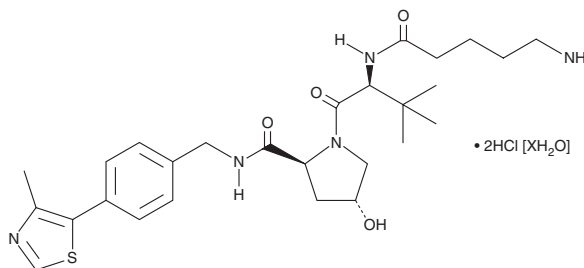
Purity: ≥95%

UV/Vis.: λ_{max}: 272 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 3 (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the VH 032 linker 3 (hydrate) in the solvent of choice, which should be purged with an inert gas. VH 032 linker 3 (hydrate) is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of VH 032 linker 3 (hydrate) in DMSO and water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

VH 032 linker 3 is a derivative of the proteolysis-targeting chimera technology (PROTAC) building block VHL ligand 1 (Item No. 21591).

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, 03/05/2021

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