

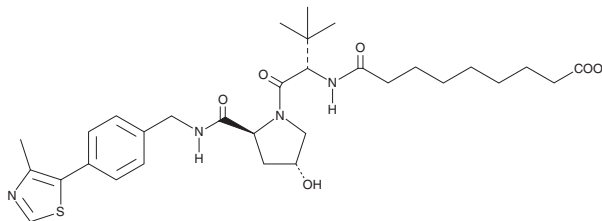
# PRODUCT INFORMATION



## VH 032 amide C7 Acid

Item No. 40447

**CAS Registry No.:** 2172819-76-8  
**Formal Name:** N-(8-carboxy-1-oxooctyl)-3-methyl-L-valyl-4R-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide  
**MF:** C<sub>31</sub>H<sub>44</sub>N<sub>4</sub>O<sub>6</sub>S  
**FW:** 600.8  
**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

VH 032 amide C7 acid is supplied as a solid. A stock solution may be made by dissolving the VH 032 amide C7 acid in the solvent of choice, which should be purged with an inert gas. VH 032 amide C7 acid is soluble in organic solvents such as ethanol and DMSO. VH 032 amide C7 acid is sparingly soluble (1-10 mg/ml) in ethanol. VH 032 amide C7 acid is soluble (≥10 mg/ml) in DMSO.

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 amide C7 acid can be prepared by directly dissolving the solid in aqueous buffers. VH 032 amide C7 acid is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

VH 032 amide C7 acid is a precursor in the synthesis of proteolysis-targeting chimeras (PROTACs) that contains a terminal carboxylic acid conjugated via a heptyl linker to VHL ligand 1 (Item No. 21591), which binds to von Hippel-Lindau (VHL) E3 ligase.<sup>1</sup> It has been used in the synthesis of PROTACs targeting anaplastic lymphoma kinase (ALK) in cancer cells.

### Reference

1. Sun, N., Ren, C., Kong, Y., *et al.* Development of a Brigatinib degrader (SIAIS117) as a potential treatment for ALK positive cancer resistance. *Eur. J. Med. Chem.* **193**, 112190 (2020).

#### 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, 09/12/2024

#### 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