

# PRODUCT INFORMATION



VH 298

Item No. 21133

Sold under license from The University of Dundee

**CAS Registry No.:** 2097381-85-4  
**Formal Name:** (2S,4R)-1-((S)-2-(1-cyanocyclopropane-1-carboxamido)-3,3-dimethylbutanoyl)-4-hydroxy-N-(4-(4-methylthiazol-5-yl)benzyl)pyrrolidine-2-carboxamide

**MF:** C<sub>27</sub>H<sub>33</sub>N<sub>5</sub>O<sub>4</sub>S

**FW:** 523.7

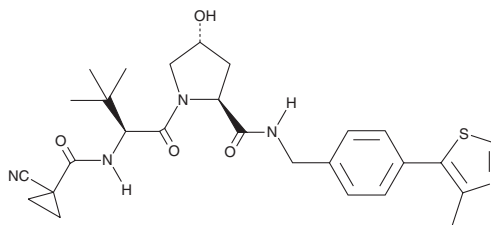
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 271 nm

**Supplied as:** A crystalline 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 298 is supplied as a crystalline solid. A stock solution may be made by dissolving the VH 298 in the solvent of choice, which should be purged with an inert gas. VH 298 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of VH 298 in these solvents is approximately 30 mg/ml.

VH 298 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VH 298 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. VH 298 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

VH 298 is a cell-permeant inhibitor of von Hippel-Lindau disease tumor suppressor (VHL; K<sub>d</sub> = 90 nM by isothermal titration calorimetry).<sup>1</sup> VHL is a component of the complex that polyubiquitinates hydroxylated hypoxia-inducible factor-α (HIF-α) isoforms leading to proteasomal degradation. Through its effects on VHL, VH 298 promotes the accumulation of HIF-α in a concentration- and time-dependent manner, resulting in the upregulation of HIF-target genes.<sup>1</sup>

## Reference

1. Frost, J.M., Galdeano, C., Soares, P., *et al.* Potent and selective chemical probe of hypoxic signalling downstream of HIF-α hydroxylation via VHL inhibition. *Nat. Commun.* **7**, 13312 (2016).

### 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, 12/15/2022

## 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