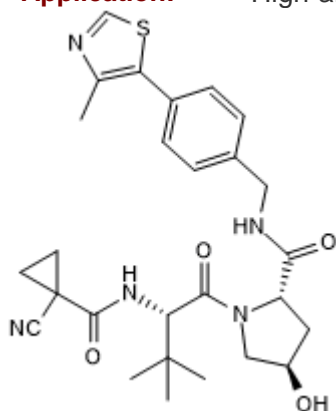


**VH298****Cat. # PC-1003****Background:**

High-affinity inhibitor of E3 ubiquitin ligase VHL ( $K_d = 80-90$  nM). Blocks interaction between VHL and HIF- $\alpha$  downstream of HIF- $\alpha$  hydroxylation, initiating hypoxic response. Results in time- and concentration-dependent accumulation of hydroxylated HIF- $\alpha$ , and upregulates mRNA and protein levels of HIF target genes. Cell permeable.

**Application:**

High-affinity inhibitor of VHL

**Product Information**

<b>Purity:</b>	>98%
<b>MW:</b>	523.65
<b>Formula:</b>	<b>C<sub>27</sub>H<sub>33</sub>N<sub>5</sub>O<sub>4</sub>S</b>
<b>CAS No.</b>	2097381-85-4
<b>Physical State:</b>	Lyophilized white powder
<b>Quantity:</b>	5mg; 10mg; 25 mg
<b>Solubility:</b>	50 mg/mL in DMSO; 50 mg/mL in Ethanol
<b>Storage:</b>	Store desiccated as supplied at -20°C for up to 3 years. Store solutions at -80°C for up to 6 months or -20°C for up to 1 month.

**References**

1. **Frost *et al*** (2015) Potent and selective chemical probe of hypoxic signalling downstream of HIF- $\alpha$  hydroxylation via VHL inhibition. *Nat. Commun.* **7** 13312 PMID: [27811928](https://pubmed.ncbi.nlm.nih.gov/27811928/)
2. **Soares *et al*** (2018) Group-based optimization of potent and cell-active inhibitors of the von Hippel-Lindau (VHL) E3 ubiquitin ligase: structure-activity relationships leading to the chemical probe (2S,4R)-1-((S)-2-(1-Cyanocyclopropanecarboxamido)-3 J. Med. Chem. **61** 599 PMID: [28853884](https://pubmed.ncbi.nlm.nih.gov/28853884/)

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2013 LifeSensors, Inc. All Rights Reserved