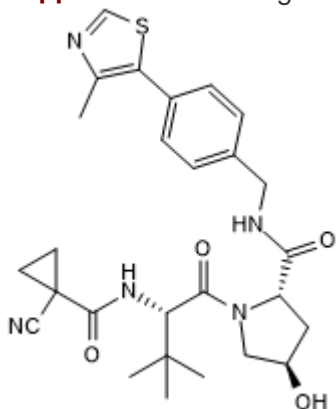


VH298**Cat. # PC-1003**

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

Application: High-affinity inhibitor of VHL

**Product Information**

Purity:	>98%
MW:	523.65
Formula:	C₂₇H₃₃N₅O₄S
CAS No.	2097381-85-4
Physical State:	Lyophilized white powder
Quantity:	5mg; 10mg; 25 mg
Solubility:	50 mg/mL in DMSO; 50 mg/mL in Ethanol
Storage:	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- α 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 (2*S*,4*R*)-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