

PRODUCT INFORMATION

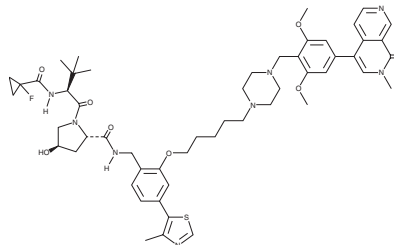


VZ 185

Item No. 41454

CAS Registry No.: 2306193-61-1
Formal Name: (2S,4R)-N-(2-((5-(4-(2,6-dimethoxy-4-(2-methyl-1-oxo-1,2-dihydro-2,7-naphthyridin-4-yl)benzyl)piperazin-1-yl)pentyl)oxy)-4-(4-methylthiazol-5-yl)benzyl)-1-((S)-2-(1-fluorocyclopropane-1-carboxamido)-3,3-dimethylbutanoyl)-4-hydroxypyrrolidine-2-carboxamide

MF: C₅₃H₆₇FN₈O₈S
FW: 995.2
Purity: ≥98%
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

VZ 185 is supplied as a solid. A stock solution may be made by dissolving the VZ 185 in the solvent of choice, which should be purged with an inert gas. VZ 185 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

VZ 185 is a proteolysis-targeting chimera (PROTAC) composed of a von Hippel-Lindau (VHL) ligand and bromodomain-containing protein 7 (BRD7)- and BRD9-binding ligand.¹ It is selective for BRD7 and BRD9 over 6,273 other proteins at 100 nM. VZ 185 degrades BRD7 and BRD9 with half-maximal degradation concentration (DC₅₀) values of 4.5 and 1.8 nM, respectively, in a cell-free assay. It decreases BRD7 and BRD9 levels in HeLa cells by 10 and 14%, respectively, when used at a concentration of 1 μM. VZ 185 reduces the proliferation of EoL-1 leukemia and A-204 sarcoma cells (IC₅₀s = 3.4 and 39 nM, respectively).

Reference

1. Zoppi, V., Hughes, S.J., Maniaci, C., *et al.* Iterative design and optimization of initially inactive proteolysis targeting chimeras (PROTACs) identify VZ185 as a potent, fast, and selective von hippel-lindau (VHL) based dual degrader probe of BRD9 and BRD7. *J. Med. Chem.* **62**(2), 699-726 (2019).

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, 02/05/2025

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