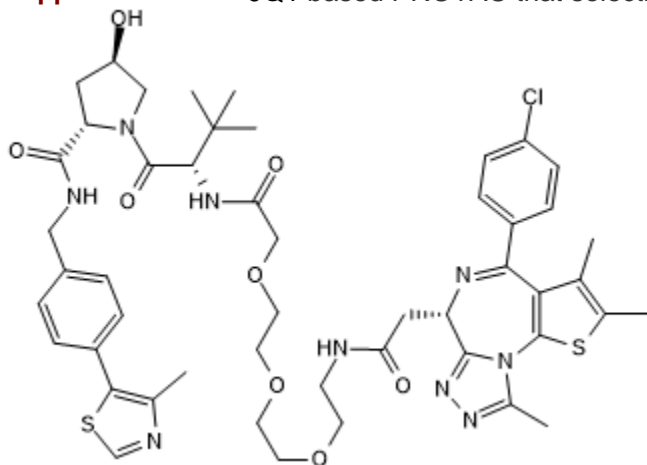


MZ1**Cat. # PC-1001****Background:**

Cell penetrant Proteolysis Targeting Chimera (PROTAC) compound based on (+)-JQ1 conjugated to a von Hippel-Lindau (VHL) ligand. Retains high affinity for BRD2, BRD3 and BRD4 bromodomains ($K_d = 13-60$ nM) but induces preferential degradation of BRD4 over BRD2 and BRD3 (DC₅₀ values for degradation of BRD4 are 8 and 23 nM in H661 and H838 cells, respectively). Exhibits potent cytotoxicity and anti-proliferative effects in AML cell lines (pEC₅₀ = 7.6 in Mv4-11 cells).

Application:

JQ1 based PROTAC that selectively degrades BRD4 in cells

**Product Information**

Purity:	>98.0%
MW:	1002.64
Formula:	C ₄₉ H ₆₀ ClN ₉ O ₈ S ₂
CAS No.	1797406-69-9
Physical State:	Lyophilized white to off-white solid
Quantity:	1mg; 5mg
Solubility:	100 mg/mL in DMSO; or 100 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. **Zengerle et al** (2015) Selective small molecule induced degradation of the BET bromodomain protein BRD4. *ACS.Chem.Biol* **10** 1770 PMID: [26035625](#)
2. **Gadd et al** (2017) Structural basis of PROTAC cooperative recognition for selective protein degradation. *Nat.Chem.Biol.* PMID: [28288108](#)
3. **Wurz et al** (2017) A "click chemistry platform" for the rapid synthesis of bispecific molecules for inducing protein degradation. *J.Med.Chem.* PMID: [28378579](#)

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses

Copyright © 2013 LifeSensors, Inc. All Rights Reserved