

UNC1999

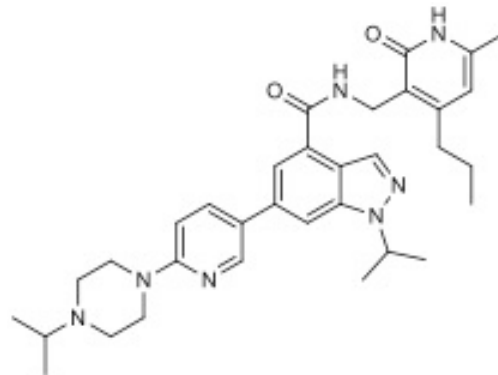
**Catalog No:** 14114, 14115**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 569.7

 $C_{33}H_{43}N_7O_2$ 

CAS 1431612-23-5

Physical Properties: White powder

**Names:** UNC1999, N[(6-methyl-2-oxo-4-propyl-1H-pyridin-3-yl)methyl]-1-propan-2-yl-6-[6-(4-propan-2-ylpiperazin-1-yl)pyridin-3-yl]indazole-4-carboxamide

**Pharmacology:** UNC1999 is an orally active, highly selective inhibitor of both wild-type and mutant EZH1 and EZH2 lysine methyltransferases ( $IC_{50}$  = 45 nM and 2 nM respectively)(ref 1). Inhibition of EZH2 with UNC1999 enhanced the efficacy of gefitinib in suppressing the proliferation of colon cancer cells (ref 2). It suppresses global H3K27 trimethylation/demethylation and inhibits the growth of mixed lineage leukemia (MLL)-rearranged leukemia cells (ref 3).

**Solubilization:** May be dissolved in DMSO (10 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**

&gt;98% (HPLC); NMR (Conforms)

## References:

1. KD Konze *et al.* *ACS Chem. Biol.* 2013, 8:1324
2. BW Katona *et al.* *Cancer Biol. Ther.* 2014, 15:1677
3. B Xu *et al.* *Blood* 2015, 125:346

**Storage and Guarantee:** Store desiccated as supplied at  $-20^{\circ}$  for up to 2 years. Store solutions at  $-20^{\circ}C$  for up to 3 months. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.