

UNC1215

Catalog #: 27405

Lot #: 121312

Size: 5 mg

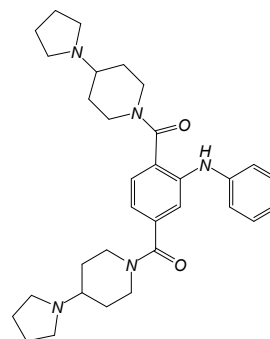
Structure:

CAS Registry #: N/A

Purity: $\geq 95\%$

Chemical Formula: $C_{32}H_{43}N_5O_2 \cdot 3TFA$

Molecular Weight: 871.8



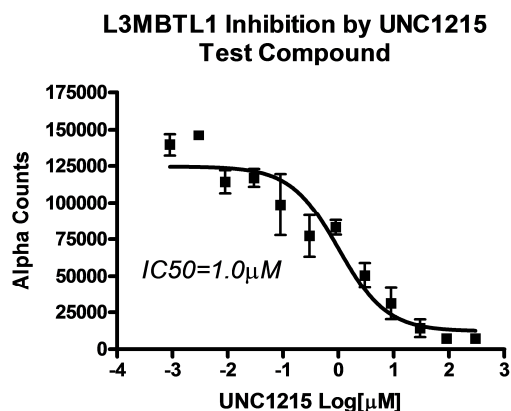
Description: UNC 1215 is a methyl-lysine inhibitor. UNC1215 inhibits L3MBTL3 with 50x more potency than other MBT family members.

Appearance: Solid.

Solubility: Soluble in water (> 10 mM). HYGROSCOPIC.

Storage/Stability: Store solid at or below -20°C ; protect from moisture. Store aqueous solutions in aliquots at -80°C for up to one month.

Biological Activity: UNC1215 has been shown to inhibit L3MBTL1 with an IC_{50} of $1.0 \mu\text{M}$.



Quality Control: The purity was determined by HPLC.

Reference: James, L.I., et al. *Nat Chem Biol.* 2013 Mar;9(3):184-91.



Data Sheet