

Product Information

Description	Tubacin is a potent, selective, reversible, and cell-permeable inhibitor of HDAC6 with an IC50 value of 4 nM.
Target:	HDAC6
Chemical Name:	N-[4-[(2R,4R,6S)-4-[(4,5-diphenyl-1,3-oxazol-2-yl)sulfanylmethyl]-6-[4-(hydroxymethyl)phenyl]-1,3-dioxan-2-yl]phenyl]-N'-hydroxyoctanediamide
Molecular Formula:	C ₄₁ H ₄₃ N ₃ O ₇ S
Supplied As:	Solid
Solubility:	insoluble in EtOH; insoluble in H ₂ O; ≥7.19 mg/mL in DMSO. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while.
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	721.86 Da
Cas No.	537049-40-4
Purity:	≥95% by HPLC

Chemical Structure

