

I-CBP112 (Hydrochloride)

Catalog #: 27336

Lot #: 140404

Size: 1 mg

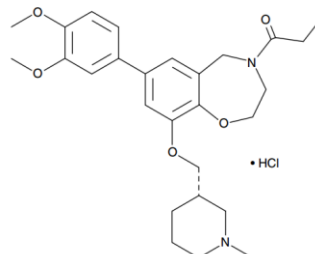
Structure:

CAS Registry #: NA

Purity: ≥98%

Chemical Formula: C₂₇H₃₆N₂O₅

Molecular Weight: 505.1

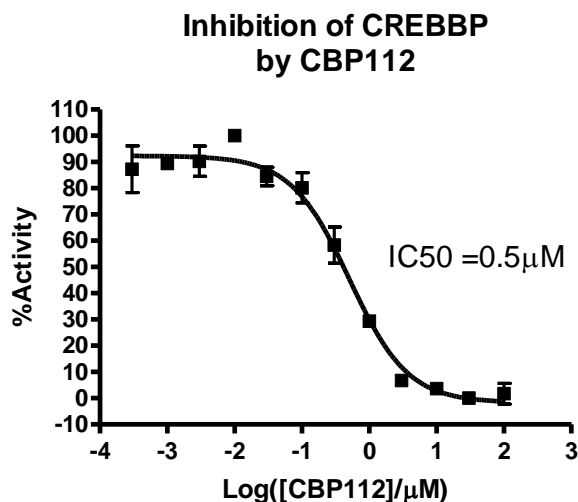


Description: I-CBP112 (Hydrochloride) is a selective inhibitor of cAMP responsive-element binding protein (CREBBP) and E1A binding protein p3000 (EP3000). The inhibitor has been shown to decrease EZH2 activity and to inhibit trimethylation of lysine 27 on histone H3 in cultured human AML HL-60 and OCI-AML3 cells. It has the same effect on primary AML cells in a dose-dependent manner.

Appearance: A crystalline solid

Solubility: Soluble in DMSO, ethanol, and dimethylformamide at approximately 16 mg/ml.

Biological Activity: I-CBP112 directly binds the bromodomains of CREBBP and EP300 with a K_D of 0.5 and 0.625 μM, respectively. It shows weak cross-reactivity with bromodomain and extra-terminal (BET) proteins and no reactivity with other bromodomain-containing proteins.





Data Sheet

Storage/Stability: Store at or below -20°C for up to two years. Protect stock solutions from moisture. Avoid freeze/thaw cycles.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Iver, N.G. *et al.*, *Oncogene*. 2004;23(**24**):4225-4231.
2. Petrij, F. *et al.*, *Nature*. 1995;376(**6538**):348-351
3. Vizmanos, J.L. *et al.*, *Gene Chromosome Canc.* 2003;**36**:402-405