

# PRODUCT INFORMATION

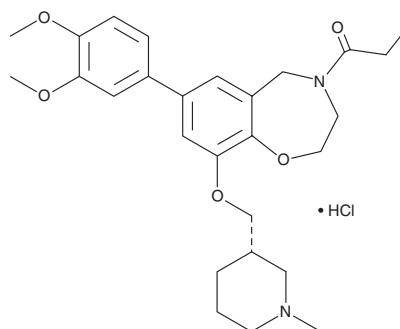


## I-CBP112 (hydrochloride)

Item No. 14468

**CAS Registry No.:** 2147701-33-3  
**Formal Name:** 1-[7-(3,4-dimethoxyphenyl)-2,3-dihydro-9-[[[(3S)-1-methyl-3-piperidiny]methoxy]-1,4-benzoxazepin-4(5H)-yl]-1-propanone, monohydrochloride

**MF:** C<sub>27</sub>H<sub>36</sub>N<sub>2</sub>O<sub>5</sub> • HCl  
**FW:** 505.1  
**Purity:** ≥90%  
**UV/Vis.:** λ<sub>max</sub>: 213, 269, 291 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

I-CBP112 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the I-CBP112 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. I-CBP112 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of I-CBP112 (hydrochloride) in these solvents is approximately 16 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of I-CBP112 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of I-CBP112 (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

I-CBP112 is an inhibitor of p300 and CREB-binding protein (CBP) histone acetyltransferases.<sup>1</sup> It binds to the p300 and CBP bromodomains (K<sub>d</sub>s = 167 and 151 nM, respectively) and is selective for p300 and CBP over BRD4, as well as a panel of 104 nuclear receptors and ion channels and a panel of 32 enzymes at 10 μM. I-CBP112 displaces acetylated histones from CBP in a cell-free assay (IC<sub>50</sub> = 170 nM). It reduces colony formation and increases differentiation of primary murine leukemic blasts and delays disease initiation following leukemic blast transplantation into sub-lethally irradiated mice when used at concentrations of 5 and 10 μM.

### Reference

1. Picaud, S., Fedorov, O., Thanasopoulou, A., *et al.* Generation of a selective small molecule inhibitor of the CBP/p300 bromodomain for leukemia therapy. *Cancer Res.* **75(23)**, 5106-5119 (2015).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/03/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM