

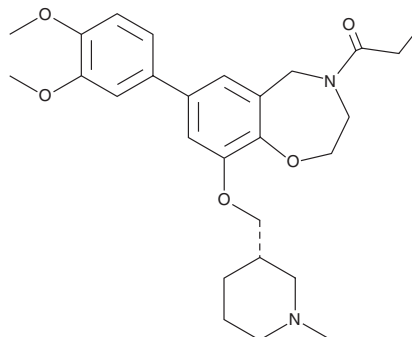
PRODUCT INFORMATION



I-CBP112

Item No. 36811

CAS Registry No.: 1640282-31-0
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
MF: C₂₇H₃₆N₂O₅
FW: 468.6
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

I-CBP112 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of I-CBP112 in these solvents is approximately 5 mg/ml in DMSO and DMF.

Description

I-CBP112 is an inhibitor of p300 and CREB-binding protein (CBP) histone acetyltransferases.¹ It binds to the p300 and CBP bromodomains (K_d 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 μ M. I-CBP112 displaces acetylated histones from CBP in a cell-free assay (IC_{50} = 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 μ M.

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, 12/21/2022

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