

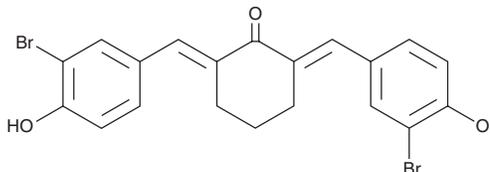
PRODUCT INFORMATION



HAT Inhibitor II

Item No. 19835

CAS Registry No.: 932749-62-7
Formal Name: 2,6-bis[(3-bromo-4-hydroxyphenyl)methylene]-cyclohexanone
Synonyms: Histone Acetyltransferase Inhibitor II
MF: C₂₀H₁₆Br₂O₃
FW: 464.2
Purity: ≥98%
UV/Vis.: λ_{max}: 249, 366 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

HAT Inhibitor II is supplied as a crystalline solid. A stock solution may be made by dissolving the HAT inhibitor II in the solvent of choice, which should be purged with an inert gas. HAT Inhibitor II is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HAT inhibitor II in these solvents is approximately 20, 30, and 50 mg/ml, respectively.

HAT Inhibitor II is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HAT inhibitor II should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HAT Inhibitor II has a solubility of approximately 0.1 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HAT Inhibitor II is a cell-permeable, selective inhibitor of the histone acetyltransferase p300 (IC₅₀ = 5 μM).¹ It does not inhibit PCAF or GCN5 acetyltransferases.¹

Reference

1. Costi, R., DiSanto, R., Artico, M., et al. Cinnamoyl compounds as simple molecules that inhibit p300 histone acetyltransferase. *J. Med. Chem.* **50(8)**, 1973-1977 (2007).

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, 10/27/2022

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