

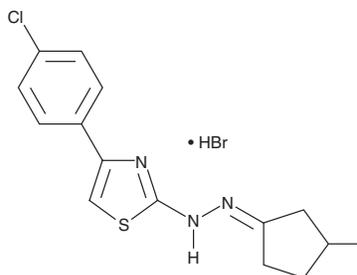
PRODUCT INFORMATION



CPTH6 (hydrobromide)

Item No. 19828

CAS Registry No.: 2321332-57-2
Formal Name: 3-methyl-cyclopentanone, 2-[4-(4-chlorophenyl)-2-thiazolyl]hydrazone, monohydrobromide
MF: C₁₅H₁₆ClN₃S • HBr
FW: 386.7
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 278 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

CPTH6 (hydrobromide) is supplied as a crystalline solid. A stock solution may be made by dissolving the CPTH6 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. CPTH6 (hydrobromide) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CPTH6 (hydrobromide) in ethanol is approximately 0.5 mg/ml and approximately 25 mg/ml in DMSO and DMF.

CPTH6 (hydrobromide) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CPTH6 (hydrobromide) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CPTH6 (hydrobromide) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CPTH6 is a thiazole derivative that inhibits the lysine acetyltransferase activity of Gcn5 and pCAF but not p300 or CBP.¹ It blocks the acetylation of H3/H4 histones and α-tubulin in several leukemia cell lines.¹ CPTH6 reduces cell viability, arresting cell cycling in the G₀/G₁ phase and inducing apoptosis.¹ It impairs autophagy in a variety of tumor cell lines, at least in part by altering ATG7-mediated elongation of autophagosomal membranes.²

References

1. Trisciuglio, D., Ragazzoni, Y., Pelosi, A., *et al.* CPTH6, a thiazole derivative, induces histone hypoacetylation and apoptosis in human leukemia cells. *Clin. Cancer Res.* **18**(2), 475-486 (2012).
2. Ragazzoni, Y., Desideri, M., Gabellini, C., *et al.* The thiazole derivative CPTH6 impairs autophagy. *Cell Death Dis.* **4**(3), e524 (2013).

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, 11/09/2022

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