

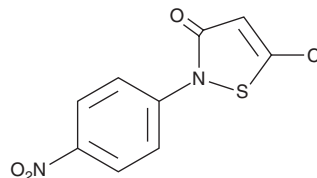
PRODUCT INFORMATION



CCT077791

Item No. 43794

CAS Registry No.: 748777-47-1
Formal Name: 5-chloro-2-(4-nitrophenyl)-3(2H)-isothiazolone
MF: C₉H₅ClN₂O₃S
FW: 256.7
Purity: ≥95%
Supplied as: A 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

CCT077791 is supplied as a solid. A stock solution may be made by dissolving the CCT077791 in the solvent of choice, which should be purged with an inert gas. CCT077791 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

CCT077791 is an inhibitor of the histone acetyltransferase p300 (IC₅₀ = 7.3 μM).¹ It is also active against *S. aureus*, *S. epidermidis*, and *B. subtilis* (MICs = 2.5, 0.3, and 2.5 μg/ml, respectively).² CCT077791 (10 μM) decreases global protein, histone H3 and H4, and α-tubulin acetylation in HCT116 colorectal cancer cells.¹ It inhibits growth of HCT116 cells and HT-29 colon cancer cells (GI₅₀s = 3 and 2.1 μM, respectively).

References

1. Stimson, L., Rowlands, M.G., Newbatt, Y.M., et al. Isothiazolones as inhibitors of PCAF and p300 histone acetyltransferase activity. *Mol. Cancer Ther.* **4(10)**, 1521-1532 (2005).
2. Khalaj, A., Adibpour, N., Shahverdi, A.R., et al. Synthesis and antibacterial activity of 2-(4-substituted phenyl)-3(2H)-isothiazolones. *Eur. J. Med. Chem.* **39(8)**, 699-705 (2004).

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, 10/08/2025

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