

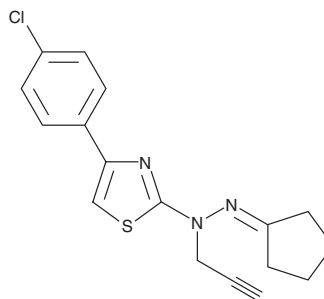
PRODUCT INFORMATION



CAY10685

Item No. 16086

CAS Registry No.: 1613116-16-7
Formal Name: cyclopentanone, 2-[4-(4-chlorophenyl)-2-thiazolyl]-2-(2-propyn-1-yl)hydrazone
Synonym: Click Tag™ CAY10685
MF: C₁₇H₁₆ClN₃S
FW: 329.8
Purity: ≥98%
UV/Vis.: λ_{max}: 249 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

CAY10685 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10685 in the solvent of choice, which should be purged with an inert gas. CAY10685 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10685 in these solvents is approximately 3, 5, and 10 mg/ml, respectively.

CAY10685 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10685 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10685 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10685 is a cell-active analog of the lysine acetyltransferase inhibitor CPTH2 (Item No. 12086) that contains an alkyne moiety for use in click chemistry reactions.¹ Its ability to inhibit N-acetyltransferase 10 activity has been used to modulate chromatin organization in order to study changes in nuclear architecture associated with cancer and certain laminopathies.¹

Reference

1. Larrieu, D., Britton, S., Demir, M., *et al.* Chemical inhibition of NAT10 corrects defects of laminopathic cells. *Science* **344**(6183), 527-532 (2014).

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, 02/06/2024

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