

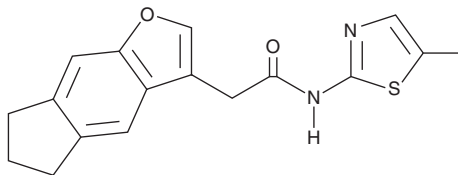
PRODUCT INFORMATION



NCT-02

Item No. 39568

CAS Registry No.: 790245-61-3
Formal Name: 6,7-dihydro-N-(5-methyl-2-thiazolyl)-5H-indeno[5,6-b]furan-3-acetamide
MF: C₁₇H₁₆N₂O₂S
FW: 312.4
Purity: ≥98%
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

NCT-02 is supplied as a solid. A stock solution may be made by dissolving the NCT-02 in the solvent of choice, which should be purged with an inert gas. NCT-02 is soluble in acetonitrile and DMSO.

Description

NCT-02 is a degrader of cyclin K.¹ It reduces the levels of cyclin K in patient-derived colorectal cancer (CRC) cells in a concentration-dependent manner. NCT-02 inhibits cancer cell proliferation in CRC tissue samples isolated from the colon (IC₅₀s = 0.9-20.6 μM), as well as from hepatic, lung, and rectal metastases (IC₅₀s = 0.5-11.8, 1, and 1.4-38.4 μM, respectively). It decreases the levels of Ataxia-telangiectasia mutated kinase (ATM) and RAD51 in patient-derived CRC tissues when used at concentrations 0.7 and 0.5 μM, respectively. NCT-02 (3 μM) reduces the number of spheroids formed and induces poly(ADP-ribose) polymerase (PARP) cleavage in patient-derived CRC spheroid cultures.

Reference

1. Dieter, S.M., Siegl, C., Codó, P.L., *et al.* Degradation of CCNK/CDK12 is a druggable vulnerability of colorectal cancer. *Cell Rep.* **36(3)**, 109394 (2021).

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, 11/17/2023

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