

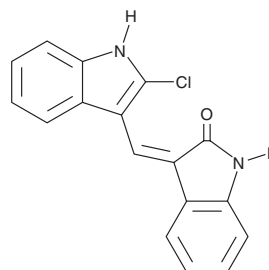
PRODUCT INFORMATION



Cdk1 Inhibitor

Item No. 21501

CAS Registry No.: 220749-41-7
Formal Name: 3-[(2-chloro-1H-indol-3-yl)methylene]-1,3-dihydro-2H-indol-2-one
Synonym: Cyclin-dependent Kinase 1 Inhibitor
MF: C₁₇H₁₁ClN₂O
FW: 294.7
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

Cdk1 inhibitor is supplied as a solid. A stock solution may be made by dissolving the Cdk1 inhibitor in the solvent of choice, which should be purged with an inert gas. Cdk1 inhibitor is soluble in the organic solvent DMSO.

Description

Cdk1 inhibitor is an inhibitor of Cdk1 (IC₅₀ = 5.8 μM).¹ It is selective for Cdk1 over Cdk5 and glycogen synthase kinase 3β (GSK3β; IC₅₀s = 25 and >100 μM, respectively). It inhibits the proliferation of HeLa cells with an IC₅₀ value of 2 μM.

Reference

1. Andreani, A., Cavalli, A., Granaiola, M., *et al.* Imidazo[2,1-b]thiazolymethylene- and indolymethylene-2-indolinones: A new class of cyclin-dependent kinase inhibitors. Design, synthesis, and CDK1/cyclin B inhibition. *Anticancer Drug Des.* **15(6)**, 447-452 (2000).

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, 12/16/2022

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