

PRODUCT INFORMATION



SKPin C1

Item No. 26141

CAS Registry No.: 432001-69-9
Formal Name: 2-[4-bromo-2-[[4-oxo-3-(3-pyridinylmethyl)-2-thioxo-5-thiazolidinylidene]methyl]phenoxy]-acetic acid

MF: C₁₈H₁₃BrN₂O₄S₂
FW: 465.3

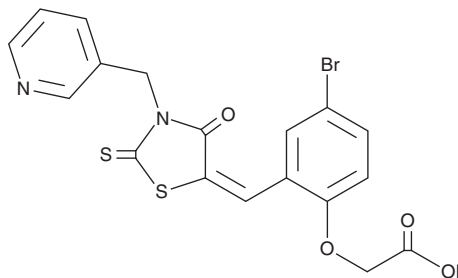
Purity: ≥98%

UV/Vis.: λ_{max}: 389 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

SKPin C1 is supplied as a crystalline solid. A stock solution may be made by dissolving the SKPin C1 in the solvent of choice. SKPin C1 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

Description

SKPin C1 is an inhibitor of Skp1-Cullin1-F-box (SCF) family protein Skp2-mediated p27 degradation.¹ It binds to Skp2 at the Cdc kinase subunit 1 (Cks1) interaction interface and blocks ubiquitylation of p27 in a Cks1-dependent manner *in vitro* when used at a concentration of 50 μM. SKPin C1 increases nuclear accumulation of p27 in ECC-1 endometrial carcinoma cells.²

References

1. Wu, L., Grigoryan, A.V., Li, Y., *et al.* Specific small molecule inhibitors of Skp2-mediated p27 degradation. *Chem. Biol.* **19**(12), 1515-1524 (2012).
2. Pavlides, S.C., Huang, K.-T., Reid, D.A., *et al.* Inhibitors of SCF-Skp2/Cks1 E3 ligase block estrogen-induced growth stimulation and degradation of nuclear p27kip1: Therapeutic potential for endometrial cancer. *Endocrinology* **154**(11), 4030-45 (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, 12/20/2022

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