

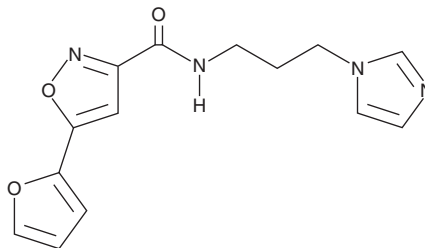
PRODUCT INFORMATION



SKL2001

Item No. 26334

CAS Registry No.: 909089-13-0
Formal Name: 5-(2-furanyl)-N-[3-(1H-imidazol-1-yl)propyl]-3-isoxazolecarboxamide
Synonym: Wnt Agonist II
MF: C₁₄H₁₄N₄O₃
FW: 286.3
Purity: ≥98%
UV/Vis.: λ_{max}: 247, 288 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

SKL2001 is supplied as a crystalline solid. A stock solution may be made by dissolving the SKL2001 in the solvent of choice. SKL2001 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of SKL2001 in these solvents is approximately 25 mg/ml.

SKL2001 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SKL2001 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SKL2001 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SKL2001 is an activator of Wnt/β-catenin signaling.^{1,2} It increases β-catenin responsive transcription in a cell-based reporter assay in a concentration-dependent manner, increases β-catenin protein levels in HEK293 cells when used at concentrations of 10 and 30 μM, and disrupts the interaction between Axin and β-catenin.¹ SKL2001 increases alkaline phosphatase (ALP) activity and mineralization in multipotent mesenchymal ST2 cells in a concentration-dependent manner, indicating increased osteoblastogenesis. SKL2001 (40 μM) reduces HCT116 and HT-29 spheroid growth and inhibits proliferation of HCT116 and HT-29 cells.² It induces cell cycle arrest at the G₀/G₁ phase and increases E-cadherin and β-catenin protein levels in HCT116 cells when used at a concentration of 40 μM.

References

1. Gwak, J., Hwang, S.G., Park, H.S., *et al.* Small molecule-based disruption of the Axin/β-catenin protein complex regulates mesenchymal stem cell differentiation. *Cell Res.* **22(1)**, 237-247 (2012).
2. Ohashi, W., Yamamine, N., Imura, J., *et al.* SKL2001 suppresses colon cancer spheroid growth through regulation of the E-cadherin/β-Catenin complex. *Biochem. Biophys. Res. Commun.* **493(3)**, 1342-1348 (2017).

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, 11/03/2022

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