

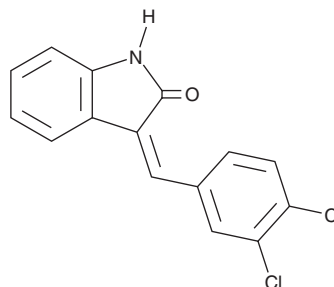
PRODUCT INFORMATION



SU 5201

Item No. 34333

CAS Registry No.: 114727-43-4
Formal Name: 3-[(3,4-dichlorophenyl)methylene]-1,3-dihydro-2H-indol-2-one
Synonym: NSC 247030
MF: C₁₅H₉Cl₂NO
FW: 290.1
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 326 nm
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

SU 5201 is supplied as a solid. A stock solution may be made by dissolving the SU 5201 in the solvent of choice, which should be purged with an inert gas. SU 5201 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SU 5201 in these solvents is approximately 20 and 30 mg/ml, respectively.

SU 5201 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SU 5201 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SU 5201 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

SU 5201 is an inhibitor of IL-2 production.¹ It inhibits IL-2 production in peripheral blood mononuclear cells (PBMCs) with an IC₅₀ value of 1 nM. SU 5201 also inhibits VEGF- or acidic FGF-induced proliferation of human umbilical vein endothelial cells (HUVECs) when used at concentrations of 2.5 and 2.3 μM, respectively.²

References

1. Bouert, E.M.L., Fensholdt, J., Nielsen, S.F., *et al.* Novel therapeutic use of indolinone derivatives. *Leo Pharma A/S. WO2005/058309AI* (2005).
2. Masferrer, J.L., Cherrington, J.M., Leahy, K.M., *et al.* Methods using a combination of a 3-heteroaryl-2-indolinone and a cyclooxygenase-2 inhibitor for the treatment of neoplasia. *Pharmacia Corporation. US2003/0119895AI* (2003).

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, 11/22/2022

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