

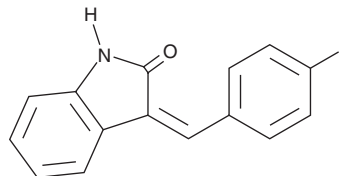
PRODUCT INFORMATION



SU 5205

Item No. 27310

CAS Registry No.: 3476-86-6
Formal Name: 3-[(4-fluorophenyl)methylene]-
1,3-dihydro-2H-indol-2-one
MF: $C_{15}H_{10}FNO$
FW: 239.2
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SU 5205 is supplied as a crystalline solid. A stock solution may be made by dissolving the SU 5205 in the solvent of choice, which should be purged with an inert gas. SU 5205 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SU 5205 in ethanol is approximately 5 mg/ml and approximately 50 mg/ml in DMSO and DMF.

SU 5205 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SU 5205 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SU 5205 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SU 5205 is an inhibitor of VEGF receptor 2 (VEGFR2/FLK1; $IC_{50} = 9.6 \mu M$).¹ It inhibits human umbilical vein endothelial cell (HUVEC) mitogenesis induced by VEGF or acidic FGF with IC_{50} values of 0.9 and 0.6 μM , respectively.

Reference

1. Tang, P.C., Sun, L., and McMahon, G. 3-(2'-halobenzylidenyl)-2-indolinone compounds for the treatment of disease. *Sugen, Inc.* **US5834504A** (1998).

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, 11/30/2022

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