

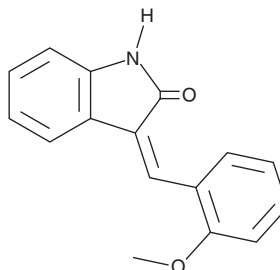
PRODUCT INFORMATION



SU 5214

Item No. 24670

CAS Registry No.: 186611-04-1
Formal Name: 1,3-dihydro-3-[(2-methoxyphenyl)methylene]-2H-indol-2-one
MF: C₁₆H₁₃NO₂
FW: 251.3
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 336 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

SU 5214 is supplied as a crystalline solid. A stock solution may be made by dissolving the SU 5214 in the solvent of choice, which should be purged with an inert gas. SU 5214 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SU 5214 in these solvents is approximately 33 mg/ml.

SU 5214 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SU 5214 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SU 5214 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 5214 is a 3-heteroaryl-2-indoline that was designed in a drug discovery program for development of tyrosine kinase inhibitors.¹

Reference

1. Tang, P.C., Sun, L., and McMahon, G. 3-heteroaryl-2-indolinone compounds for the treatment of disease. *Sugen, Inc. US005792783A* (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, 04/20/2023

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