

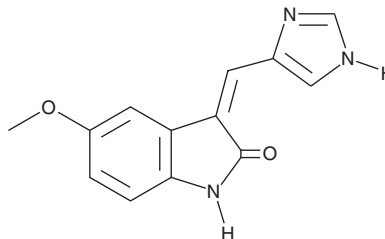
# PRODUCT INFORMATION



## SU 9516

Item No. 14796

**CAS Registry No.:** 377090-84-1  
**Formal Name:** 1,3-dihydro-3Z-(1H-imidazol-5-ylmethylene)-5-methoxy-2H-indol-2-one  
**MF:** C<sub>13</sub>H<sub>11</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 241.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 271, 291, 348, 354 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 9516 is supplied as a crystalline solid. A stock solution may be made by dissolving the SU 9516 in the solvent of choice, which should be purged with an inert gas. SU 9516 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SU 9516 in these solvents is approximately 5 mg/ml.

SU 9516 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SU 9516 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SU 9516 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

1. Lane, M.E., Yu, B., Rice, A., *et al.* A novel cdk2-selective inhibitor, SU9516, induces apoptosis in colon carcinoma cells. *Cancer Res.* **61(16)**, 6170-6177 (2001).
2. Yu, B., Lane, M.E., and Wadler, S. SU9516, a cyclin-dependent kinase 2 inhibitor, promotes accumulation of high molecular weight E2F complexes in human colon carcinoma cells. *Biochem. Pharmacol.* **64(7)**, 1091-1100 (2002).
3. Gao, N., Kramer, L., Rahmani, M., *et al.* The three-substituted indolinone cyclin-dependent kinase 2 inhibitor 3-[1-(3H-imidazol-4-yl)-meth-(Z)-ylidene]-5-methoxy-1,3-dihydro-indol-2-one (SU9516) kills human leukemia cells *via* down-regulation of Mcl-1 through a transcriptional mechanism. *Mol. Pharmacol.* **70(2)**, 645-655 (2006).

#### 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, 11/09/2022

#### 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