

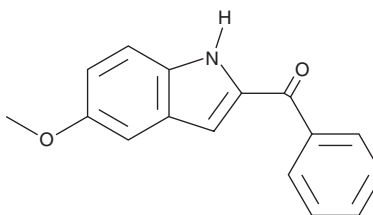
PRODUCT INFORMATION



D-64131

Item No. 22944

CAS Registry No.: 74588-78-6
Formal Name: (5-methoxy-1H-indol-2-yl)phenylmethanone
MF: C₁₆H₁₃NO₂
FW: 251.3
Purity: ≥98%
UV/Vis.: λ_{max}: 324 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

D-64131 is supplied as a crystalline solid. A stock solution may be made by dissolving the D-64131 in the solvent of choice, which should be purged with an inert gas. D-64131 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of D-64131 in these solvents is approximately 25 and 33 mg/ml, respectively.

D-64131 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, D-64131 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. D-64131 has a solubility of approximately 0.16 mg/ml in a 1:10 solution of DMF:PBS (pH 7.0) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

D-64131 is an antimitotic agent that inhibits polymerization of microtubules with an IC₅₀ value of 0.53 μM in a radioligand binding assay using purified tubules from bovine brain.¹ It arrests the cell cycle at the G₂/M transition (IC₅₀ = 92 nM) and is preferentially cytotoxic to cycling cells. It has antiproliferative activity in cancer cell lines derived from various tissues and organs (IC₅₀s = 24-144 nM) and in cell lines resistant to other tubulin inhibitors (IC₅₀s = 20-77 nM). D-64131 (200 and 400 mg/kg) is also efficacious in reducing tumor growth in a mouse xenograft model of human amelanotic melanoma.

Reference

1. Beckers, T., Reissmann, T., Schmidt, M., *et al.* 2-aryloxyindoles, a novel class of potent, orally active small molecule tubulin inhibitors. *Cancer Res.* **62(11)**, 3113-3119 (2002).

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, 11/07/2022

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