

PRODUCT INFORMATION



Stauprimide

Item No. 15398

CAS Registry No.: 154589-96-5
Formal Name: N-[(9S,10R,11R,13R)-2,3,10,11,12,13-hexahydro-10-methoxy-9-methyl-1,3-dioxo-9,13-epoxy-1H,9H-diindolo[1,2,3-gh:3',2',1'-lm]pyrrolo[3,4-j][1,7]benzodiazonin-11-yl]-N-methylbenzamide

Synonym: N-Benzoyl-7-oxo Staurosporine

MF: C₃₅H₂₈N₄O₅

FW: 584.6

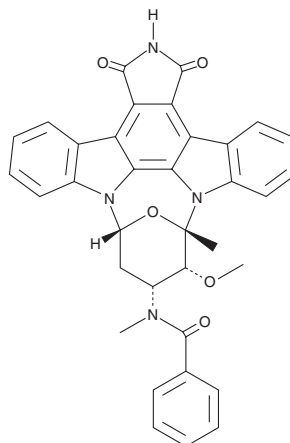
Purity: ≥98%

UV/Vis.: λ_{max}: 239, 287, 317, 412 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

Stauprimide is supplied as a crystalline solid. A stock solution may be made by dissolving the stauprimide in the solvent of choice. Stauprimide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of stauprimide in these solvents is approximately 0.2, 10, and 25 mg/ml, respectively.

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

Stauprimide is a small molecule that increases the efficiency of mouse and human embryonic stem cell differentiation in conjunction with lineage specification cues (EC₅₀ = 0.3 μM).¹ It specifically inhibits the nuclear localization of NME2, which results in the suppression of c-Myc, a key regulator of pluripotency.¹

Reference

1. Zhu, S., Wurdak, H., Wang, J., *et al.* A small molecule primes embryonic stem cells for differentiation. *Cell Stem Cell* 4(5), 416-426 (2009).

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, 11/16/2022

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