

PRODUCT INFORMATION



1,2,3,4-Tetrahydrostaurosporine

Item No. 29987

CAS Registry No.: 220038-19-7
Formal Name: (9S,10R,11R,13R)-2,3,10,11,12,13,15,16,17,18-decahydro-10-methoxy-9-methyl-11-(methylamino)-9,13-epoxy-1H,9H-diindolo[1,2,3-gh:3',2',1'-lm]pyrrolo[3,4-j][1,7]benzodiazonin-1-one

Synonym: AFN-941

MF: C₂₈H₃₀N₄O₃

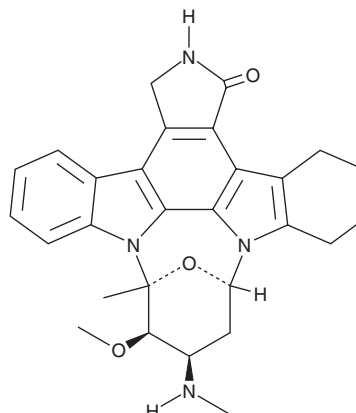
FW: 470.6

Purity: ≥95%

Supplied as: A 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

1,2,3,4-Tetrahydrostaurosporine is supplied as a solid. A stock solution may be made by dissolving the 1,2,3,4-tetrahydrostaurosporine in the solvent of choice, which should be purged with an inert gas. 1,2,3,4-Tetrahydrostaurosporine is slightly soluble in acetone and DMSO.

Description

1,2,3,4-Tetrahydrostaurosporine is a derivative of staurosporine (Item No. 81590) and an inhibitor of mutant EGFR (IC₅₀ = 74 nM for EGFR^{T790M}).¹ It is selective for EGFR^{T790M} over wild-type EGFR (IC₅₀ = 390 nM). It also binds to Janus kinase 3 (JAK3).²

References

1. Song, X., Liu, X., and Ding, X. Staurosporine scaffold-based rational discovery of the wild-type sparing reversible inhibitors of EGFR T790M gatekeeper mutant in lung cancer with analog-sensitive kinase technology. *J. Mol. Recognit.* **30(4):e2590** (2017).
2. Boggon, T.J., Li, Y., Manley, P.W., et al. Crystal structure of the Jak3 kinase domain in complex with a staurosporine analog. *Blood* **106(3)**, 996-1002 (2005).

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, 12/08/2022

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