

PRODUCT INFORMATION

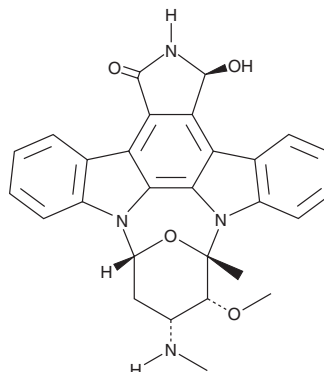


UCN-02

Item No. 20221

CAS Registry No.: 121569-61-7
Formal Name: (3S,9S,10R,11R,13R)-2,3,10,11,12,13-hexahydro-3-hydroxy-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: 7-*epi*-hydroxy Staurosporine

MF: C₂₈H₂₆N₄O₄
FW: 482.5
Purity: ≥95%
Supplied as: A 1 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

UCN-02 is a derivative of staurosporine (Item No. 81590) and a stereoisomer of UCN-01 (Item No. 18130) first isolated from a high staurosporine-producing *Streptomyces* culture as a minor co-metabolite.¹ UCN-02 inhibits protein kinase C with a slightly reduced potency than UCN-01 (IC₅₀ = 62 and 4.1 nM, for UCN-02 and UCN-01, respectively) and is cytotoxic to the growth of HeLa S3 cells.¹

Reference

1. Takahashi, I., Saitoh, Y., Yoshida, M., *et al.* UCN-01 and UCN-02, new selective inhibitors of protein kinase C. II. Purification, physico-chemical properties, structural determination and biological activities. *J. Antibiot.* **42**(4), 571-576 (1989).

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, 12/15/2025

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