

PRODUCT INFORMATION



Rubratoxin A

Item No. 19605

CAS Registry No.: 22467-31-8
Formal Name: (4S,5R,10R)-*rel*-10-[(R)-[(2R)-3,6-dihydro-6-oxo-2H-pyran-2-yl]hydroxymethyl]-4,5,8,9,10,11-hexahydro-4,8-dihydroxy-5-[[1S]-1-hydroxyheptyl]-1H-cyclonona[1,2-c:5,6-c']difuran-1,3,6-trione

Synonyms: BRN 1633008, HSDB 3532

MF: C₂₆H₃₂O₁₁

FW: 520.5

Purity: ≥95%

Supplied as: A powder

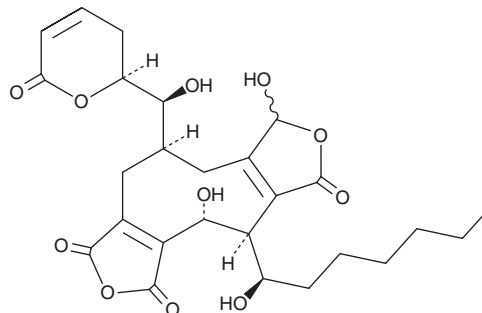
Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Penicillium* sp.

Special Conditions: Keep cool and dry

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



Laboratory Procedures

Rubratoxin A is supplied as a powder. A stock solution may be made by dissolving the rubratoxin A in the solvent of choice, which should be purged with an inert gas. Rubratoxin A is soluble in acetonitrile and DMSO. Rubratoxin A is unstable in alcohols, such as methanol and ethanol, and may cause esterification.

Description

Rubratoxin A is a natural mycotoxin and competitive inhibitor of protein phosphatase 2A (PP2A; IC₅₀ = 170 nM).¹ It has no significant effect on several other phosphatases, including PP1, PP2B, and protein tyrosine phosphatase 1B. Rubratoxin A drives overphosphorylation of PP2A substrate proteins in cultured cells treated for three hours at 20 μM.¹ It causes suppression of tumor metastasis and reduction of primary tumor volume in mouse xenografts.¹

Reference

1. Wada, S., Usami, I., Umezawa, Y., *et al.* Rubratoxin A specifically and potently inhibits protein phosphatase 2A and suppresses cancer metastasis. *Cancer Sci.* **101**(3), 743-750 (2010).

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, 08/09/2023

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