

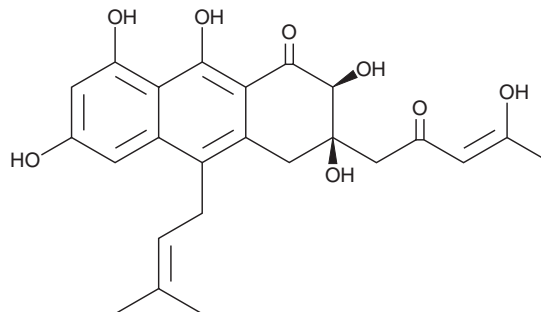
PRODUCT INFORMATION



Neosartoricin B

Item No. 33162

CAS Registry No.: 1421708-43-1
Formal Name: (2S,3R)-3,4-dihydro-2,3,6,8,9-pentahydroxy-3-[(3Z)-4-hydroxy-2-oxo-3-penten-1-yl]-10-(3-methyl-2-buten-1-yl)-1(2H)-anthracenone
Synonym: NSC 34758
MF: C₂₄H₂₆O₈
FW: 442.5
Purity: ≥85%
UV/Vis.: λ_{max}: 234, 278 nm
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

Neosartoricin B is supplied as a solid. A stock solution may be made by dissolving the neosartoricin B in the solvent of choice, which should be purged with an inert gas. Neosartoricin B is soluble in DMSO.

Description

Neosartoricin B is a polyketide produced in *A. nidulans* heterologously expressing dermatophyte-derived polyketide synthase genes.¹

Reference

1. Yin, W.-B., Chooi, Y.H., Smith, A.R., et al. Discovery of cryptic polyketide metabolites from dermatophytes using heterologous expression in *Aspergillus nidulans*. *ACS Synth. Biol.* **2**(11), 629-634 (2013).

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, 11/17/2022

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