

# PRODUCT INFORMATION

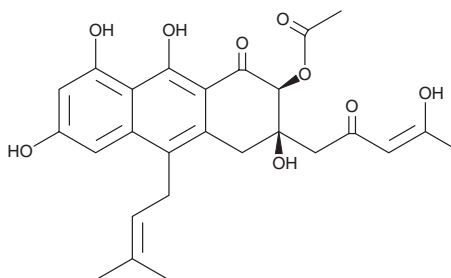


## Neosartoricin

Item No. 33103

**CAS Registry No.:** 1421941-29-8  
**Formal Name:** (2S,3R)-2-(acetyloxy)-3,4-dihydro-3,6,8,9-tetrahydro-3-[(3Z)-4-hydroxy-2-oxo-3-penten-1-yl]-10-(3-methyl-2-buten-1-yl)-1(2H)-anthracenone

**MF:** C<sub>26</sub>H<sub>28</sub>O<sub>9</sub>  
**FW:** 484.5  
**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

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

### Description

Neosartoricin is a prenylated tricyclic polyketide produced in *A. fumigatus* and *N. fischeri* overexpressing the *N. fischeri* polycyclic prenyltransferase (pcPTase) *nsc* gene cluster.<sup>1</sup> It inhibits the proliferation of anti-CD3 and anti-CD28-stimulated mouse spleen T cells (IC<sub>50</sub> = 2.99 μM).

### Reference

1. Chooi, Y.-H., Fang, J., Liu, H., *et al.* Genome mining of a prenylated and immunosuppressive polyketide from pathogenic fungi. *Org. Lett.* **15**(4), 780-783 (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, 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