

PRODUCT INFORMATION



16-*epi* Latrunculin B

Item No. 19608

CAS Registry No.: 444911-05-1
Formal Name: (4S)-4-[(1R,4Z,8Z,10S,13R,15R)-15-hydroxy-5,10-dimethyl-3-oxo-2,14-dioxabicyclo[11.3.1]heptadeca-4,8-dien-15-yl]-2-thiazolidinone

MF: C₂₀H₂₉NO₅S

FW: 395.5

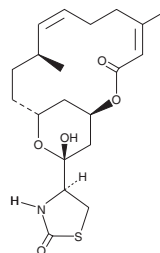
Purity: ≥95%

Supplied as: A film

Storage: -20°C

Stability: ≥4 years

Item Origin: Sponge/*Negombata magnifica*



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

Laboratory Procedures

16-*epi* Latrunculin B is supplied as a solid. A stock solution may be made by dissolving the 16-*epi* latrunculin B in the solvent of choice. 16-*epi* Latrunculin B is soluble in organic solvents such as ethanol and DMSO.

Description

16-*epi* Latrunculin B, first isolated from the Red Sea sponge *N. magnifica*, is a stereoisomer of the actin polymerization inhibitor, latrunculin B (Item No. 10010631).^{1,2} At 5-10 µg/ml, 16-*epi* latrunculin B can disrupt microfilament activity in an actin disruption assay.² It also demonstrates cytotoxic activity in mouse KA31T and NIH3T3 tumor cells (GI₅₀S = 1 and 4 µg/ml, respectively) and antiviral activity against herpes simplex type 1 virus (ED₅₀ = 1 µg/ml).¹

References

- Hoye, T.R., Ayyad, S.-E.N., Eklov, B.M., *et al.* Toward computing relative configurations: 16-*epi*-Latrunculin B, a new stereoisomer of the actin polymerization inhibitor latrunculin B. *J. Am. Chem. Soc.* **124**(25), 7405-7410 (2016).
- Vilozny, B., Amagata, T., Mooberry, S.L., *et al.* A new dimension to the biosynthetic products isolated from the sponge *Negombata magnifica*. *J. Nat. Prod.* **67**(6), 1055-1057 (2004).

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, 10/26/2022

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