

PRODUCT INFORMATION



Bafilomycin D

Item No. 19438

CAS Registry No.: 98813-13-9

Formal Name: (3Z,5E,7R,8S,9S,11E,13E,15S,16R)-16-
[(1S,2R,3S,5E,7S,8R)-2,8-dihydroxy-
1,3,7,9-tetramethyl-4-oxo-5-decen-1-
yl]-8-hydroxy-3,15-dimethoxy-5,7,9,11-
tetramethyl-oxacyclohexadeca-3,5,11,13-
tetraen-2-one

MF: C₃₅H₅₆O₈

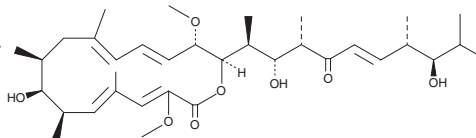
FW: 604.8

Purity: ≥99%

Supplied as: A powder

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Bafilomycin D is supplied as a powder. A stock solution may be made by dissolving the bafilomycin D in the solvent of choice, which should be purged with an inert gas. Bafilomycin D is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Bafilomycin D is a bacterial metabolite that has been found in *Streptomyces* and has diverse biological activities.¹⁻³ It is an inhibitor of vacuolar H⁺-ATPases (V-ATPases; K_i = 20 nM in *N. crassa* vacuolar membranes) and is selective for V-ATPases over P-type ATPases (K_i = 20,000 nM for the *E. coli* enzyme).² Bafilomycin D induces autophagosome accumulation in MCF-7 cells when used at concentrations ranging from 10 to 1,000 nM.¹ It reduces viral genome copy numbers in the culture supernatant of Vero E6 cells infected with H1N1 influenza A.³

References

1. Carr, G., Williams, D.E., Díaz-Marrero, A.R., *et al.* Bafilomycins produced in culture by *Streptomyces* spp. isolated from marine habitats are potent inhibitors of autophagy. *J. Nat. Prod.* **73**(3), 422-427 (2010).
2. Dröse, S., Bindseil, K.U., Bowman, E.J., *et al.* Inhibitory effect of modified bafilomycins and concanamycins on P- and V-type adenosinetriphosphatases. *Biochemistry* **32**(15), 3902-3906 (1993).
3. Xie, X., Lu, S., Pan, X., *et al.* Antiviral bafilomycins from a feces-inhabiting *Streptomyces* sp. *J. Nat. Prod.* **84**(2), 537-543 (2021).

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, 01/26/2024

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