

PRODUCT INFORMATION

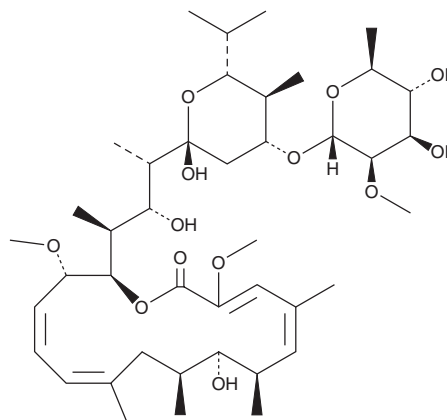


Leucanicidin

Item No. 25479

CAS Registry No.: 91021-66-8
Formal Name: (3Z,5E,7R,8S,9S,11E,13E,15S,16R)-16-[[[(1S,2R,3S)-3-[(2R,4R,5S,6R)-4-[(6-deoxy-2-O-methyl- α -L-mannopyranosyl)oxy]tetrahydro-2-hydroxy-5-methyl-6-(1-methylethyl)-2H-pyran-2-yl]-2-hydroxy-1-methylbutyl]-8-hydroxy-3,15-dimethoxy-5,7,9,11-tetramethyl-oxacyclohexadeca-3,5,11,13-tetraen-2-one

MF: C₄₂H₇₀O₁₃
FW: 783.0
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Leucanicidin is supplied as a solid. A stock solution may be made by dissolving the leucanicidin in the solvent of choice, which should be purged with an inert gas. Leucanicidin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. NOTE: Forms methyl ketal on long term storage in methanol.

Leucanicidin is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Leucanicidin is a macrolide bacterial metabolite originally isolated from *S. halstedii*.¹ It is toxic to *L. separata* fourth instar larvae when used at a concentration of 20 ppm and to *H. contortus*, *T. colubriformis*, and *O. circumcincta* larvae (LD₅₀s = 0.23-0.42 μ g/ml).²

References

1. Isogai, A., Sakuda, S., Matsumoto, S., et al. The structure of leucanicidin, a novel insecticidal macrolide produced by *Streptomyces halstedii*. *Agr. Biol. Chem.* **48(5)**, 1379-1381 (1984).
2. Lacey, E., Gill, J.H., Power, M.L., et al. Bafilolides, potent inhibitors of the motility and development of the free-living stages of parasitic nematodes. *Int. J. Parasitol.* **25(3)**, 349-357 (1995).

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, 11/28/2022

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