

PRODUCT INFORMATION



IB-96212

Item No. 40938

CAS Registry No.: 220858-11-7
Formal Name: (4E,18E,20E)-3',4',5',6'-tetrahydro-7,9,11,12,13,14,15-heptahydroxy-6'-(2-hydroxy-1-methylbutyl)-22-(2-hydroxypropyl)-5',6,8,16,28,29-hexamethyl-14-propyl-10-[[tetrahydro-5-hydroxy-6-methyl-2H-pyran-2-yl]oxy]-spiro[2,26-dioxabicyclo[23.3.1]nonacosan-4,18,20-triene-27,2'-[2H]pyran]-3-one

MF: C₅₄H₉₄O₁₆

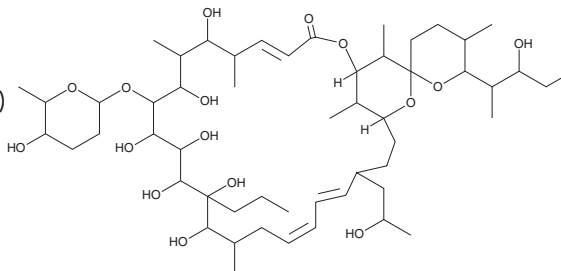
FW: 999.3

Purity: ≥70%

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

IB-96212 is supplied as a solid. A stock solution may be made by dissolving the IB-96212 in the solvent of choice, which should be purged with an inert gas. IB-96212 is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

IB-96212 is a macrolide that has been found in *Micromonospora* and has antibacterial and anticancer activities.¹ It is active against *M. luteus*, *B. subtilis*, and *S. aureus* (MICs = 0.4, 100, and 100 µg/ml, respectively). IB-96212 induces death in P388 leukemia, A549 lung cancer, HT-29 colon cancer, and SK-MEL-28 melanoma cells (IC₅₀s = 0.0001, 1, 1, and 1 µg/ml, respectively).

Reference

1. Fernández-Chimeno, R.I., Cañedo, L., Espliego, F., *et al.* IB-96212, a novel cytotoxic macrolide produced by a marine *Micromonospora*. I. Taxonomy, fermentation, isolation and biological activities. *J. Antibiot. (Tokyo)* **53**(5), 474-478 (2000).

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, 07/02/2024

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