

PRODUCT INFORMATION



18-Deoxyherboxidiene

Item No. 29063

CAS Registry No.: 1200128-66-0
Formal Name: (2R,5S,6S)-tetrahydro-6-[(1E,3E,5S)-6-[(2R,3R)-3-[(1R,2S)-2-methoxy-1-methylbutyl]-2-methyl-2-oxiranyl]-1,5-dimethyl-1,3-hexadien-1-yl]-5-methyl-2H-pyran-2-acetic acid

Synonym: RQN-18690A

MF: C₂₅H₄₂O₅

FW: 422.6

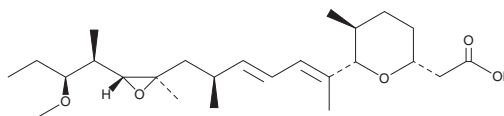
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

18-Deoxyherboxidiene is supplied as a solid. A stock solution may be made by dissolving the 18-deoxyherboxidiene in the solvent of choice, which should be purged with an inert gas. 18-Deoxyherboxidiene is soluble in ethanol, methanol, and DMSO.

Description

18-Deoxyherboxidiene is a bacterial metabolite that has been found in *Streptomyces* QN18690 and has antiangiogenic activity.¹ It binds to the SF3b subunit of spliceosome-associated protein 130 (SAP130), SAP145, or SAP155 in HeLa cell extracts when used at 100, 1,000, and 10,000 ng and induces the accumulation of unspliced mRNA in HEK293T cells in a reporter assay. 18-Deoxyherboxidiene inhibits VEGF-induced migration of human umbilical vein endothelial cells (HUVECs) in a concentration-dependent manner and VEGF-induced HUVEC tube formation when used at a concentration of 0.3 µg/ml.

Reference

1. Kakeya, H., Kaida, D., Sekiya, H., *et al.* RQN-18690A (18-deoxyherboxidiene) targets SF3b, a spliceosome component, and inhibits angiogenesis. *J. Antibiot. (Tokyo)* **69(2)**, 121-123 (2016).

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, 10/17/2022

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