

PRODUCT INFORMATION



13-acetyl-9-Dihydrobaccatin III

Item No. 29701

CAS Registry No.: 142203-65-4
Formal Name: (2aR,4S,4aS,5R,6R,9S,11S,12S,12aR,12bS)-2a,3,4,4a,5,6,9,10,12,12a-decahydro-4a,8,13,13-tetramethyl-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxete-4,5,6,9,11,12,12b-heptol

Synonym: 9-Dihydro-13-acetyl-baccatin III

MF: C₃₃H₄₂O₁₂

FW: 630.7

Purity: ≥95%

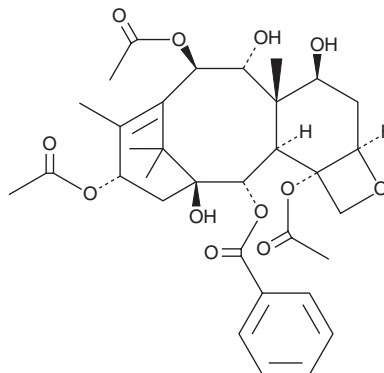
UV/Vis.: λ_{max}: 229 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Taxus chinensis*



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

Laboratory Procedures

13-acetyl-9-Dihydrobaccatin III is supplied as a solid. A stock solution may be made by dissolving the 13-acetyl-9-dihydrobaccatin III in the solvent of choice, which should be purged with an inert gas. 13-acetyl-9-Dihydrobaccatin III is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of 13-acetyl-9-dihydrobaccatin III in these solvents is approximately 5 and 10 mg/ml, respectively.

13-acetyl-9-Dihydrobaccatin III is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 13-acetyl-9-dihydrobaccatin III should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 13-acetyl-9-Dihydrobaccatin III has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

13-acetyl-9-Dihydrobaccatin III is a taxane originally isolated from *T. canadensis*.¹ It inhibits proliferation of P388 leukemia cells with an IC₅₀ value of 20 µg/ml.

Reference

1. Gunawardana, G.P., Premachandran, U., Burren, N.S., et al. Isolation of 9-Dihydro-13-acetyl-baccatin III from *Taxus canadensis*. *J. Nat. Prod.* **55**(11), 1686-1689 (1992).

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, 12/08/2022

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