

PRODUCT INFORMATION



10-acetyl Docetaxel

Item No. 34497

CAS Registry No.: 125354-16-7
Formal Name: (αR,βS)-β-[[[(1,1-dimethylethoxy)carbonyl]amino]-α-hydroxy-benzenepropanoic acid, (2aR,4S,4aS,6R,9S,11S,12S,12aR,12bS)-6,12b-bis(acetyloxy)-12-(benzoyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-4,11-dihydroxy-4a,8,13,13-tetramethyl-5-oxo-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester PNU 101383, 10-acetyl Taxotere

Synonyms:

MF: C₄₅H₅₅NO₁₅

FW: 849.9

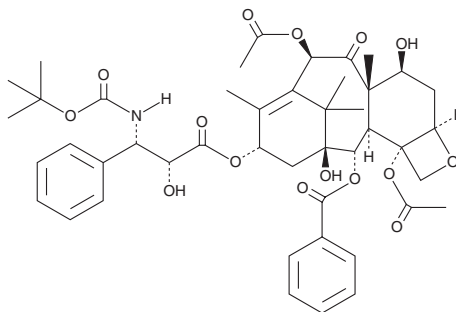
Purity: ≥98%

UV/Vis.: λ_{max}: 229 nm

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

10-acetyl Docetaxel is supplied as a solid. A stock solution may be made by dissolving the 10-acetyl docetaxel in the solvent of choice, which should be purged with an inert gas. 10-acetyl Docetaxel is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 10-acetyl docetaxel in these solvents is approximately 10, 3, and 2 mg/ml, respectively.

Description

10-acetyl Docetaxel is a derivative of paclitaxel (Item No. 10461) and an inhibitor of microtubule depolymerization.¹ It inhibits calcium-induced depolymerization of calf brain microtubule protein (MTP) in a cell-free assay when used at 20 μM. 10-acetyl Docetaxel is cytotoxic to J774.2 cells (EC₅₀ = 0.05 μM).

Reference

1. Swindell, C.S., Heering, J.M., Krauss, N.E., *et al.* Characterization of the taxol structure-activity profile for the locus of the A-ring side chain. *Bioorg. Med. Chem. Lett.* **4**(12), 1531-1536 (1994).

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, 11/22/2022

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