

PRODUCT INFORMATION



6 α -hydroxy Paclitaxel

Item No. 10009027

CAS Registry No.: 153212-75-0
Formal Name: β S-(benzoylamino)- α R-hydroxy-(2aR,3S,4R,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-3,4,11-trihydroxy-4a,8,13,13-tetramethyl-5-oxo-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester-benzenepropanoic acid

Synonym: 6 α -hydroxy Taxol

MF: C₄₇H₅₁NO₁₅

FW: 869.9

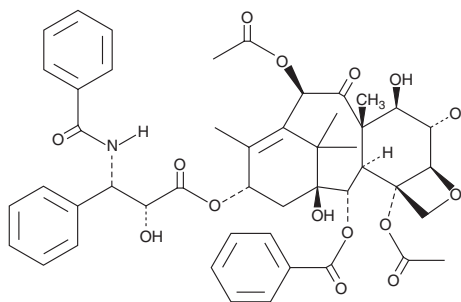
Purity: \geq 95%

UV/Vis.: λ_{\max} : 228 nm

Supplied as: A crystalline 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

6 α -hydroxy Paclitaxel is supplied as a crystalline solid. A stock solution may be made by dissolving the 6 α -hydroxy paclitaxel in the solvent of choice, which should be purged with an inert gas. 6 α -hydroxy Paclitaxel is soluble in organic solvents such as methanol and DMSO. The solubility of 6 α -hydroxy paclitaxel in these solvents is approximately 10 and 30 mg/ml, respectively.

Description

6 α -hydroxy Paclitaxel is a metabolite of the antimetabolic and anticancer agent paclitaxel (Item No. 10461).¹ It is formed from paclitaxel by the cytochrome P450 (CYP) isoform CYP2C8.^{1,2}

References

1. Kumar, G.N., Oatis, J.E., Jr., Thornburg, K.R., *et al.* 6 α -Hydroxytaxol: Isolation and identification of the major metabolite of taxol in human liver microsomes. *Drug Metab. Dispos.* **22**, 177-179 (1994).
2. Walle, T., Kumar, G.N., McMillian, J.M., *et al.* Taxol metabolism in rat hepatocytes. *Biochem. Pharmacol.* **46(9)**, 1661-1664 (1993).

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, 02/28/2023

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