

PRODUCT INFORMATION

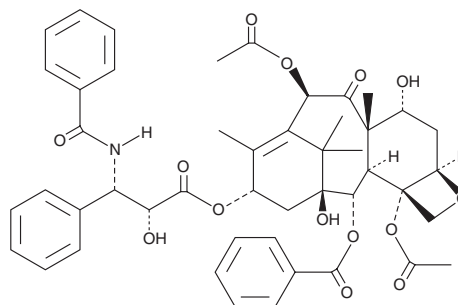


7-*epi* Paclitaxel

Item No. 20741

CAS Registry No.: 105454-04-4
Formal Name: (αR,βS)-β-(benzoylamino)-α-hydroxy-benzenepropanoic acid, (2aR,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-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

Synonym: 7-*epi* Taxol
MF: C₄₇H₅₁NO₁₄
FW: 853.9
Purity: ≥98%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A crystalline 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

7-*epi* Paclitaxel is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-*epi* paclitaxel in the solvent of choice, which should be purged with an inert gas. 7-*epi* Paclitaxel is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 7-*epi* paclitaxel in these solvents is approximately 5, 10, and 25 mg/ml, respectively.

7-*epi* Paclitaxel is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 7-*epi* paclitaxel should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 7-*epi* Paclitaxel has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

7-*epi* Paclitaxel is an active metabolite of paclitaxel (Item No. 10461).¹ It is formed from paclitaxel by hydrolysis. 7-*epi* Paclitaxel inhibits microtubule polymerization in J774.2 and CHO cells (EC₅₀ = 120 and 310 nM, respectively).²

References

1. Royer, I., Alvinerie, P., Armand, J.P., et al. Paclitaxel metabolites in human plasma and urine: Identification of 6 alpha-hydroxytaxol, 7-epitaxol and taxol hydrolysis products using liquid chromatography/atmospheric-pressure chemical ionization mass spectrometry. *Rapid Commun. Mass Spectrom.* **9**(6), 495-502 (1995).
2. Ringel, I. and Horwitz, S.B. Taxol is converted to 7-epitaxol, a biologically active isomer, in cell culture medium. *J. Pharmacol. Exp. Ther.* **242**(2), 692-698 (1987).

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, 09/14/2023

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