

PRODUCT INFORMATION



7-*epi*-10-Deacetyltaxol

Item No. 34296

CAS Registry No.: 78454-17-8

Formal Name: (α R, β S)- β -(benzoylamino)- α -hydroxy-benzenepropanoic acid, (2aR,4R,4aS,6R,9S,11S,12S,12aR,12bS)-12b-(acetyloxy)-12-(benzoyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-4,6,11-trihydroxy-4a,8,13,13-tetramethyl-5-oxo-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester

Synonyms: 10-Deacetyl-7-epitaxol, 10-Deacetyl-7-*epi*-paclitaxel, 10-Deacetyl-7-epipaclitaxel, 10-Desacetyl-7-*epi*-taxol, 10-Desacetyl-7-epipaclitaxel, 7-*epi*-10-deacetyl Paclitaxel, Ormosin VI

MF: $C_{45}H_{49}NO_{13}$

FW: 811.9

Purity: $\geq 95\%$

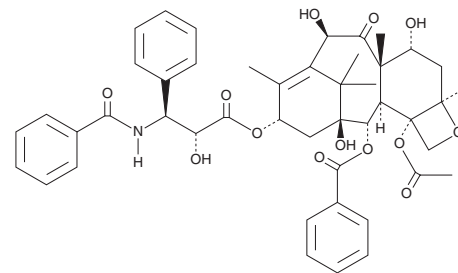
UV/Vis.: λ_{max} : 227 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Synthetic



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

Laboratory Procedures

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

Description

7-*epi*-10-Deacetyltaxol is a taxane diterpenoid and derivative of paclitaxel (Item No. 10461) that has been found in *Taxus* and has α -glucosidase inhibitory and anticancer activities.^{1,2} It inhibits α -glucosidase ($IC_{50} = 48.8 \mu\text{M}$) and is cytotoxic to HeLa cells ($IC_{50} = 0.085 \text{ nM}$).¹ 7-*epi*-10-Deacetyltaxol induces apoptosis and DNA fragmentation in HepG2 hepatocellular carcinoma cells in a concentration-dependent manner.²

References

1. Dang, P.H., Nguyen, H.X., Duong, T.T.T., *et al.* α -Glucosidase inhibitory and cytotoxic taxane diterpenoids from the stem bark of *Taxus wallichiana*. *J. Nat. Prod.* **80**(4), 1087-1095 (2017).
2. Subban, K., Singh, S., Subramani, R., *et al.* Fungal 7-*epi*-10-deacetyltaxol produced by an endophytic *Pestalotiopsis microspora* induces apoptosis in human hepatocellular carcinoma cell line (HepG2). *BMC Complement. Altern. Med.* **17**(1), 504 (2017).

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, 11/08/2022

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