

# PRODUCT INFORMATION



## 10-deacetyl Paclitaxel

Item No. 28503

**CAS Registry No.:** 78432-77-6  
**Formal Name:** ( $\alpha$ R, $\beta$ S)- $\beta$ -(benzoylamino)- $\alpha$ -hydroxy-benzenepropanoic acid, (2aR,4S,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-DAP, 10-Deacetyltaxol

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

**FW:** 811.9

**Purity:**  $\geq 98\%$

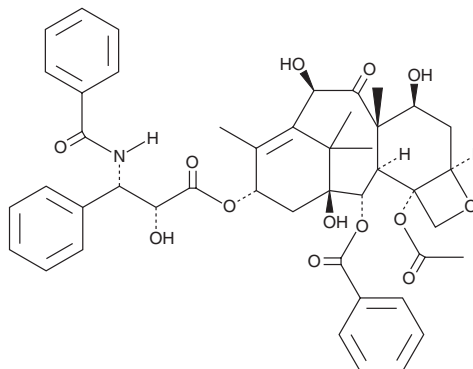
**UV/Vis.:**  $\lambda_{max}$ : 227 nm

**Supplied as:** A solid

**Storage:**  $-20^{\circ}\text{C}$

**Stability:**  $\geq 4$  years

**Item Origin:** Plant/*Taxus chinensis* (Pilger) Rehd



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

### Laboratory Procedures

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 10-deacetyl paclitaxel can be prepared by directly dissolving the solid in aqueous buffers. 10-deacetyl Paclitaxel is slightly soluble in PBS, pH 7.2. We do not recommend storing the aqueous solution for more than one day.

### Description

10-deacetyl Paclitaxel (10-DAP) is a taxane and precursor in the semisynthetic synthesis of paclitaxel (Item No. 10461) that has been found in *Taxus* and has anticancer activity.<sup>1,2</sup> It is cytotoxic to KB carcinoma cells ( $EC_{50} = 0.027 \mu\text{g/ml}$ ).<sup>1</sup> 10-DAP (16 mg/kg) reduces tumor growth in the PS mouse lymphocytic leukemia model.

### References

- McLaughlin, J.L., Miller, R.W., Powell, R.G., *et al.* 19-Hydroxybaccatin III, 10-deacetylcephalomannine, and 10-deacetyltaxol: New antitumor taxanes from *Taxus wallichiana*. *J. Nat. Prod.* **44**(3), 312-319 (1981).
- Pyo, S.H., Choi, H.J., and Han, B.H. Large-scale purification of 13-dehydroxybaccatin III and 10-deacetylpaclitaxel, semi-synthetic precursors of paclitaxel, from cell cultures of *Taxus chinensis*. *J. Chromatogr. A.* **1123**(1), 15-21 (2006).

#### 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, 12/16/2022

#### 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](http://WWW.CAYMANCHEM.COM)