

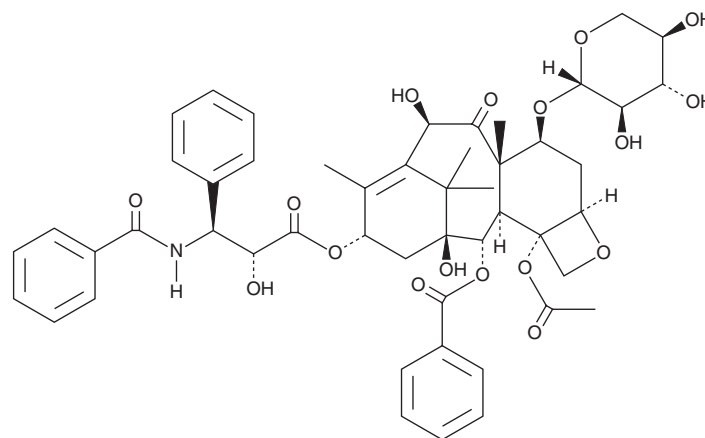
PRODUCT INFORMATION



10-deacetyl-7-xylosyl Paclitaxel

Item No. 32955

CAS Registry No.: 90332-63-1
Formal Name: (α R, β S)- β -(benzoylamino)- α -hydroxybenzenepropanoic acid, (2aR,4S,4aS,6R,9S,11S,12S,12aR,12bS)-12b-(acetyloxy)-12-(benzyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-6,11-dihydroxy-4a,8,13,13-tetramethyl-5-oxo-4-(β -D-xylopyranosyloxy)-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester
MF: C₅₀H₅₇NO₁₇
FW: 944.0
Purity: \geq 60%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

10-deacetyl-7-xylosyl Paclitaxel is supplied as a solid. A stock solution may be made by dissolving the 10-deacetyl-7-xylosyl paclitaxel in the solvent of choice, which should be purged with an inert gas. 10-deacetyl-7-xylosyl Paclitaxel is slightly soluble in DMSO (sonicated).

Description

10-deacetyl-7-xylosyl Paclitaxel is a disruptor of microtubules and derivative of paclitaxel (Item No. 10461) that has been found in *T. chinensis*.^{1,2} It induces the disassembly of pig brain and *Physarum* microtubules in cell-free assays (IC₅₀s = 0.3 and 0.5 μ M, respectively).¹ 10-deacetyl-7-xylosyl Paclitaxel inhibits proliferation of (IC₅₀ = 5 μ M), as well as induces cell cycle arrest at the G₂/M phase and apoptosis in, PC3 prostate cancer cells.²

References

1. Lataste, H., Senilh, V., Wright, M., *et al.* Relationships between the structures of taxol and baccatine III derivatives and their *in vitro* action on the disassembly of mammalian brain and *Physarum* amoebal microtubules. *Proc. Natl. Acad. Sci. USA* **81(13)**, 4090-4094 (1984).
2. Jiang, S., Zu, Y., Fu, Y., *et al.* Activation of the mitochondria-driven pathway of apoptosis in human PC-3 prostate cancer cells by a novel hydrophilic paclitaxel derivative, 7-xylosyl-10-deacetylpaclitaxel. *Int. J. Oncol.* **33(1)**, 103-111 (2008).

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, 11/10/2022

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