

PRODUCT INFORMATION



7-Xylosyltaxol

Item No. 32956

CAS Registry No.: 90332-66-4
Formal Name: (αR,βS)-β-(benzoylamino)-α-hydroxy-benzenepropanoic acid, (2aR,4S,4aS,6R,9S,11S,12,12aR,12bS)-6,12b-bis(acetyloxy)-12-(benzoyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-11-hydroxy-4a,8,13,13-tetramethyl-5-oxo-4-(D-xylopyranosyloxy)-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester

Synonym: 7-Xylosylpaclitaxel

MF: C₅₂H₅₉NO₁₈

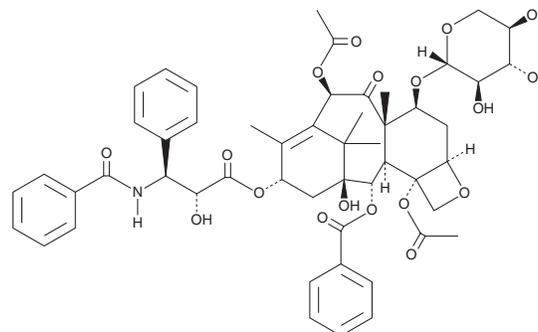
FW: 986.0

Purity: ≥98%

Supplied as: A 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-Xylosyltaxol is supplied as a solid. A stock solution may be made by dissolving the 7-xylosyltaxol in the solvent of choice, which should be purged with an inert gas. 7-Xylosyltaxol is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

7-Xylosyltaxol is a taxane and derivative of paclitaxel (Item No. 10461) that has been found in *T. chinensis* and has microtubule disruptor and anticancer activities.¹⁻³ It induces the disassembly of pig brain and *Physarum* microtubules in cell-free assays (IC₅₀s = 0.2 and 0.4 μM, respectively).² 7-Xylosyltaxol inhibits proliferation of A549, MCF-7, A2780, HCT-8, and SW480 cancer cells with IC₅₀ values of 0.39, 0.16, 0.85, 5.9, and 5.6 μM, respectively.³

References

1. Yu, S.-S., Sun, Q.-W., Zhang, X.-P., *et al.* Content and distribution of active components in cultivated and wild *Taxus chinensis* var. *mairei* plants. *Ying Yong Sheng Tai Xue Bao* **23(10)**, 2641-2647 (2012).
2. 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. U.S.A.* **81(13)**, 4090-4094 (1984).
3. Sun, Z.-H., Chen, Y., Guo, Y.-Q., *et al.* Isolation and cytotoxicity evaluation of taxanes from the barks of *Taxus wallichiana* var. *mairei*. *Bioorg. Med. Chem. Lett.* **25(6)**, 1240-1243 (2015).

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