

PRODUCT INFORMATION



Paclitaxel C

Item No. 34905

CAS Registry No.: 153415-45-3

Formal Name: (αR,βS)-α-hydroxy-β-[(1-oxohexyl)amino]-benzenepropanoic acid, (2aR,4S,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

Synonyms: Taxol C, Taxuyunnanine A

MF: C₄₆H₅₇NO₁₄

FW: 847.9

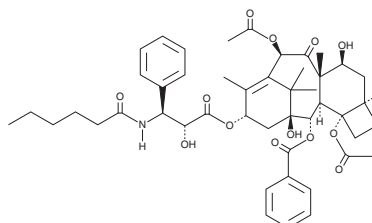
Purity: ≥95%

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

Paclitaxel C is supplied as a solid. A stock solution may be made by dissolving the paclitaxel C in the solvent of choice, which should be purged with an inert gas. Paclitaxel C is soluble in methanol.

Description

Paclitaxel C is a taxoid and derivative of paclitaxel (Item No. 10461) that has been found in *T. cuspidata* and an inhibitor of microtubule depolymerization.¹ It inhibits calcium-induced microtubule depolymerization *in vitro* when used at a concentration of 10 μM.

References

1. Kobayashi, J., Hosoyama, H., Wang, X.-X., *et al.* Effects of taxoids from *Taxus cuspidata* on microtubule depolymerization and vincristine accumulation in MDR cells. *Bioorg. Med. Chem. Lett.* **7**(4), 393-398 (1997).

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, 12/08/2022

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