

PRODUCT INFORMATION



7-acetyl Paclitaxel

Item No. 39288

CAS Registry No.: 92950-39-5

Formal Name: (αR,βS)-β-(benzoylamino)-α-hydroxy-benzenepropanoic acid, (2aR,4S,4aS,6R,9S,11S,12S,12aR,12bS)-4,6,12b-tris(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-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester

Synonym: 7-acetyl Taxol

MF: C₄₉H₅₃NO₁₅

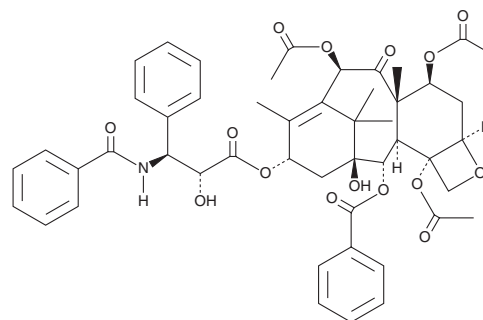
FW: 895.9

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

Description

7-acetyl Paclitaxel is an inhibitor of microtubule depolymerization.¹ It inhibits calcium-induced microtubule depolymerization when used at a concentration of 10 μM. 7-acetyl Paclitaxel reduces the growth of J774.2 macrophages (IC₅₀ = ~60 nM). It also increases the levels of nitric oxide (NO) and TNF in isolated mouse peritoneal macrophages.²

References

1. Mellado, W., Magri, N.F., Kingston, D.G., *et al.* Preparation and biological activity of taxol acetates. *Biochem. Biophys. Res. Commun.* **124**(2), 329-336 (1984).
2. Ojima, I., Fumero-Oderda, C.L., Kuduk, S.D., *et al.* Structure-activity relationship study of taxoids for their ability to activate murine macrophages as well as inhibit the growth of macrophage-like cells. *Bioorg. Med. Chem.* **11**(13), 2867-2888 (2003).

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, 09/11/2024

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