

PRODUCT INFORMATION



10-oxo Docetaxel

Item No. 21672

CAS Registry No.: 167074-97-7

Formal Name: (αR,βS)-β-[[[(1,1-dimethylethoxy)carbonyl]amino]-α-hydroxy-benzenepropanoic acid, (2aR,4S,4aS,9S,11S,12S,12aR,12bS)-12b-(acetyloxy)-12-(benzoyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-4,11-dihydroxy-4a,8,13,13-tetramethyl-5,6-dioxo-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester

MF: C₄₃H₅₁NO₁₄

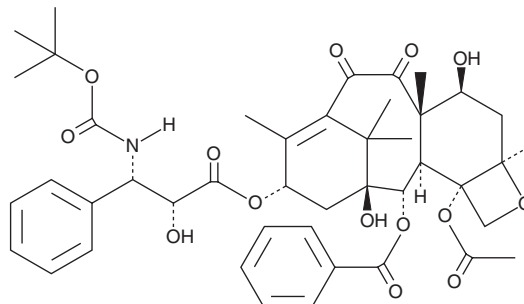
FW: 805.9

Purity: ≥85%

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

10-oxo Docetaxel is supplied as a solid. A stock solution may be made by dissolving the 10-oxo docetaxel in the solvent of choice, which should be purged with an inert gas. 10-oxo Docetaxel is slightly soluble in chloroform and methanol.

Description

10-oxo Docetaxel is a degradation product of docetaxel (Item No. 11637).¹

Reference

1. Malleswara Reddy, A., Banda, N., Govind Dagdu, S., *et al.* Evaluation of the pharmaceutical quality of docetaxel injection using new stability indicating chromatographic methods for assay and impurities. *Sci. Pharm.* **78(2)**, 215-231 (2010).

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