

PRODUCT INFORMATION



Paclitaxel octadecanedioate

Item No. 28635

CAS Registry No.: 2089211-45-8
Formal Name: octadecanedioic acid-1-[(1R,2S)-2-(benzoylamino)-1-[[[(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]oxy]carbonyl]-2-phenylethyl] ester

Synonyms: 1,18-Octadecanedioic Acid-Paclitaxel, ODDA-PTX, PTX-FA18

MF: C₆₅H₈₃NO₁₇

FW: 1,150.4

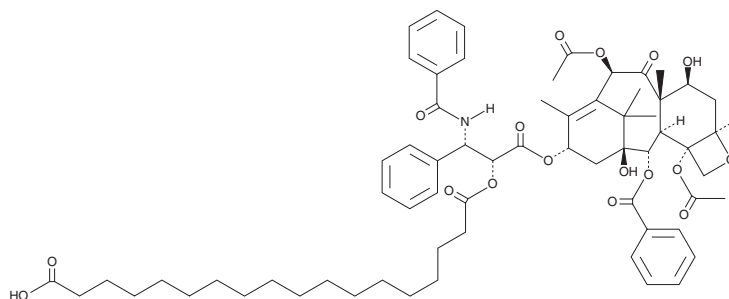
Purity: ≥98%

UV/Vis.: λ_{max}: 228 nm

Supplied as: A crystalline 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

Paclitaxel octadecanedioate is supplied as a crystalline solid. A stock solution may be made by dissolving the paclitaxel octadecanedioate in the solvent of choice, which should be purged with an inert gas. Paclitaxel octadecanedioate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of paclitaxel octadecanedioate in these solvents is approximately 20, 25, and 30 mg/ml, respectively.

Paclitaxel octadecanedioate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, paclitaxel octadecanedioate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Paclitaxel octadecanedioate has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Paclitaxel octadecanedioate is a prodrug form of paclitaxel (Item No. 10461) that is comprised of paclitaxel conjugated to 1,18-octadecanedioic acid.¹ Unlike paclitaxel, it does not promote tubulin polymerization *in vitro* when used at a concentration of 10 μM. A 5:1 mixture of paclitaxel octadecanedioate:human serum albumin (HSA) is cytotoxic to HT-1080, PANC-1, and HT-29 cells (IC₅₀s = 12, 2.48, and 8.62 nM, respectively). This mixture reduces tumor growth and increases survival in an HT-1080 mouse xenograft model in a dose-dependent manner.

Reference

1. Callmann, C.E., Leguyader, C.L.M., Burton, S.T., *et al.* Antitumor activity of 1,18-octadecanedioic acid-paclitaxel complexed with human serum albumin. *J. Am. Chem. Soc.* **141**(30), 11765-11769 (2019).

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, 03/19/2024

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