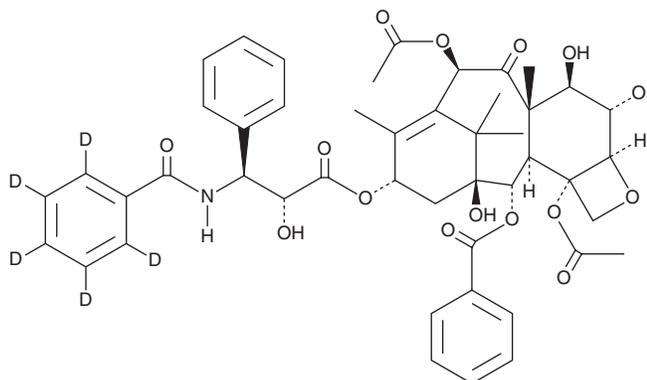


PRODUCT INFORMATION



6 α -hydroxy Paclitaxel-d₅ Item No. 40926

CAS Registry No.: 1315376-90-9
Formal Name: (α R, β S)- β -(benzoyl-2,3,4,5,6-d₅-amino)- α -hydroxy-benzenepropanoic acid, (2aR,3S,4R,4aS,6R,9S,11S,12S,12aR,12bS)-6,12b-bis(acetyloxy)-12-(benzyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-3,4,11-trihydroxy-4a,8,13,13-tetramethyl-5-oxo-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester
Synonym: 6 α -hydroxy Taxol-d₅
MF: C₄₇H₄₆D₅NO₁₅
FW: 874.9
Chemical Purity: \geq 98% (6 α -hydroxy Paclitaxel)



Deuterium Incorporation: \geq 99% deuterated forms (d₁-d₅); \leq 1% d₀
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

6 α -hydroxy Paclitaxel-d₅ is intended for use as an internal standard for the quantification of 6 α -hydroxy Paclitaxel (Item No. 10009027) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

6 α -hydroxy Paclitaxel-d₅ is supplied as a solid. A stock solution may be made by dissolving the 6 α -hydroxy Paclitaxel-d₅ in the solvent of choice, which should be purged with an inert gas. 6 α -hydroxy Paclitaxel-d₅ is soluble in methanol, DMSO, and dimethyl formamide.

Description

6 α -hydroxy Paclitaxel is a metabolite of the antimetabolic and anticancer agent paclitaxel (Item No. 10461).¹ It is formed from paclitaxel by the cytochrome P450 (CYP) isoform CYP2C8.^{1,2}

References

1. Kumar, G.N., Oatis, J.E., Jr., Thornburg, K.R., *et al.* 6 alpha-Hydroxytaxol: Isolation and identification of the major metabolite of taxol in human liver microsomes. *Drug Metab. Dispos.* **22**, 177-179 (1994).
2. Walle, T., Kumar, G.N., McMillian, J.M., *et al.* Taxol metabolism in rat hepatocytes. *Biochem. Pharmacol.* **46(9)**, 1661-1664 (1993).

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, 04/26/2024

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