

PRODUCT INFORMATION

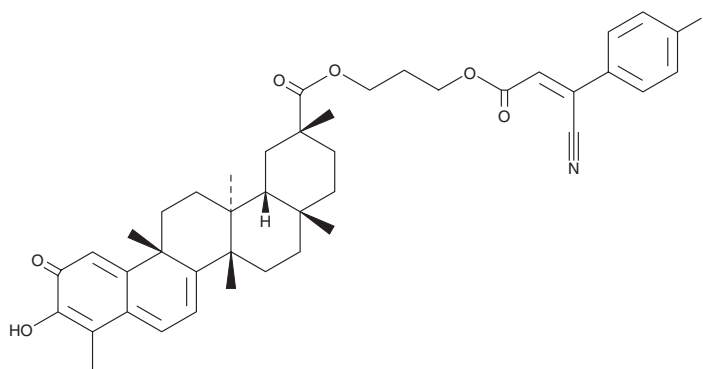


CAY10747

Item No. 29915

CAS Registry No.: 2413849-81-5
Formal Name: (9 β ,13 α ,14 β ,20 α)-3-hydroxy-9,13-dimethyl-2-oxo-24,25,26-trinoroleana-1(10),3,5,7-tetraen-29-oic acid, 3-[[[(2Z)-3-cyano-3-(4-fluorophenyl)-1-oxo-2-propen-1-yl]oxy]propyl ester

MF: C₄₂H₄₈FNO₆
FW: 681.8
Purity: \geq 98%
UV/Vis.: λ_{max} : 298, 422 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

CAY10747 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10747 in the solvent of choice, which should be purged with an inert gas. CAY10747 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10747 in these solvents is approximately 1, 20, and 30 mg/ml, respectively.

Description

CAY10747 is an inhibitor of the protein-protein interaction between heat shock protein 90 (Hsp90) and cell division cycle 37 (Cdc37) and a derivative of celastrol (Item No. 70950).¹ It decreases protein levels of the Hsp90-Cdc37 complex and the Hsp90-Cdc37 clients phosphorylated Akt and Cdk4 in A549 cells when used at a concentration of 5 μ M. CAY10747 inhibits proliferation of A549, MCF-7, HOS, and HepG2 cells (IC₅₀s = 0.41, 0.64, 0.9, and 0.94 μ M, respectively) and induces apoptosis in A549 cells.

Reference

1. Li, N., Xu, M., Wang, B., *et al.* Discovery of novel celastrol derivatives as Hsp90-Cdc37 interaction disruptors with antitumor activity. *J. Med. Chem.* **62**(23), 10798-10815 (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, 02/21/2024

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