

# PRODUCT INFORMATION

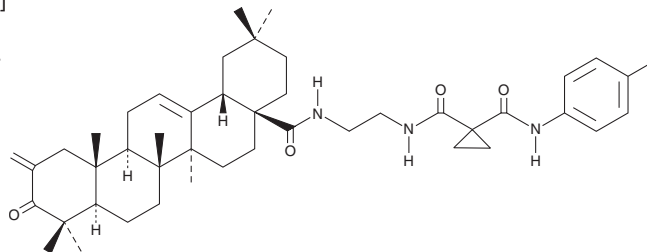


**CAY10735**

Item No. 26753

**CAS Registry No.:** 2251753-58-7  
**Formal Name:** N-[2-[[[1-[[[4-fluorophenyl]amino]carbonyl]cyclopropyl]carbonyl]amino]ethyl]-2-methylene-3-oxoolean-12-en-28-amide

**MF:** C<sub>44</sub>H<sub>60</sub>FN<sub>3</sub>O<sub>4</sub>  
**FW:** 714.0  
**Purity:** ≥90%  
**UV/Vis.:** λ<sub>max</sub>: 243 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

CAY10735 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10735 in the solvent of choice, which should be purged with an inert gas. CAY10735 is soluble in chloroform and ethanol.

## Description

CAY10735 is an anticancer compound.<sup>1</sup> It inhibits proliferation in a panel of melanoma and breast, pancreatic, and lung cancer cell lines (IC<sub>50</sub>s = 0.674-11.56 μM). CAY10735 inhibits migration of and the epithelial-to-mesenchymal transition (EMT) in A375 and B16/F10 melanoma cells *in vitro* in a concentration-dependent manner. It reduces the viability of spheroid A375 and B16/F10 cells (IC<sub>50</sub>s = 3.04 and 1.24 μM, respectively) and increases production of reactive oxygen species (ROS) in these cells in a concentration-dependent manner. CAY10735 (5 mg/kg) reduces tumor growth in B16/F10 melanoma and Lewis lung carcinoma mouse models and an A375 mouse xenograft model.

## Reference

1. Liu, X., Li, B., Zhang, Z., *et al.* Synthesis and discovery novel anti-cancer stem cells compounds derived from the natural triterpenoid acids. *J. Med. Chem.* **61**(23), 10814-10833 (2018).

### 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, 02/16/2024

## 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