

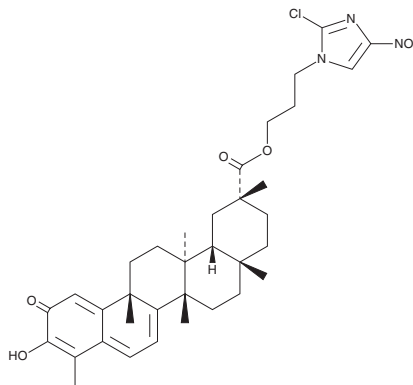
PRODUCT INFORMATION



CAY10792

Item No. 37353

CAS Registry No.: 2361009-68-7
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-(2-chloro-4-nitro-1H-imidazol-1-yl)propyl ester
MF: C₃₅H₄₄ClN₃O₆
FW: 638.2
Purity: $\geq$ 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: $\geq$ 2 years



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

Laboratory Procedures

CAY10792 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10792 in the solvent of choice, which should be purged with an inert gas. CAY10792 is soluble in chloroform.

Description

CAY10792 is an anticancer agent.¹ It binds to heat shock protein 90 (Hsp90) and cell division cycle 37 (Cdc37) in cell-free assays. CAY10792 inhibits the growth of A549, HCT116, U2OS, and MDA-MB-231 cancer cells (IC₅₀s = 0.54, 0.59, 0.57, and 0.57 μ M, respectively) but not A549 cells overexpressing Hsp90, Cdc37, or Hsp90 and Cdc37 (IC₅₀s = >100 μ M for all). It induces cell cycle arrest at the G₀/G₁ phase and apoptosis in A549 lung cancer cells when used at a concentration of 0.8 μ M. *In vivo*, CAY10792 (0.5 and 1 mg/kg) reduces tumor weight in an A549 mouse xenograft model.

Reference

1. Li, N., Xu, M., Zhang, L., *et al.* Discovery of novel celastrol-imidazole derivatives with anticancer activity *in vitro* and *in vivo*. *J. Med. Chem.* **65**(6), 4578-4589 (2022).

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, 11/11/2022

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