

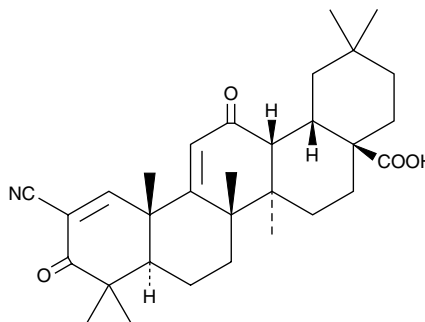
Product Information



CDDO

Item No. 81035

CAS Registry No.: 218600-44-3
Formal Name: 2-cyano-3,12-dioxo-oleana-1,9(11)-dien-28-oic acid
Synonyms: Bardoxolone, RTA 401
MF: C₃₁H₄₁NO₄
FW: 491.7
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 240 nm



Laboratory Procedures

For long term storage, we suggest that CDDO be stored as supplied at -20°C. It should be stable for at least two years.

CDDO is supplied as a crystalline solid. A stock solution may be made by dissolving the CDDO in the solvent of choice. CDDO is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of CDDO in these solvents is approximately 5 mg/ml.

CDDO is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CDDO should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CDDO has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

CDDO is a synthetic oleanane triterpenoid that blocks the cellular synthesis of inducible nitric oxide synthase and inducible COX-2 in INF-γ-activated mouse macrophages with an IC₅₀ value of 0.4 nM.¹ By suppressing reactive oxygen and nitrogen species (ROS/RNS) formation, it promotes the cellular control of ROS/RNS levels that would lead to DNA damage associated with tumorigenesis.² In various cancer cell lines, CDDO has been shown to specifically inhibit proliferation and induce apoptosis.² Mechanism studies revealed that CDDO is a ligand for peroxisome proliferator-activated receptor γ, and also that it induces genes regulated by Nrf2, including heme oxygenase-1 and eotaxin-1, which play a role in antioxidant response element signaling activity.³⁻⁵

References

1. Honda, T., Rounds, B.V., Gribble, G.W., *et al.* Design and synthesis of 2-cyano-3,12-dioxoolean-1,9-dien-28-oic acid, a novel and highly active inhibitor of nitric oxide production in mouse macrophages. *Bioorg. Med. Chem. Lett.* **8**, 2711-2714 (1998).
2. Suh, N., Wang, Y., Honda, T., *et al.* A novel synthetic oleanane triterpenoid, 2-cyano-3,12-dioxoolean-1,9-dien-28-oic acid, with potent differentiating, antiproliferative, and anti-inflammatory activity. *Cancer Res.* **59**, 336-341 (1999).
3. Ferguson, H.E., Kulkarni, A., Lehmann, G.M., *et al.* Electrophilic peroxisome proliferator-activated receptor-γ ligands have potent antifibrotic effects in human lung fibroblasts. *Am. J. Respir. Cell Mol. Biol.* **41**, 722-730 (2009).
4. Liby, K., Hock, T., Yore, M.M., *et al.* The synthetic triterpenoids, CDDO and CDDO-imidazolidine, are potent inducers of heme oxygenase-1 and Nrf2/ARE signaling. *Cancer Res.* **65**, 4789-4798 (2005).
5. Fourtounis, J., Wang, I.-M., Mathieu, M.-C., *et al.* Gene expression profiling following NRF2 and KEAP1 siRNA knockdown in human lung fibroblasts identifies CCL11/Eotaxin-1 as a novel NRF2 regulated gene. *Respir. Res.* **13**(1), [In press] (2012).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/81035

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 10/30/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com