

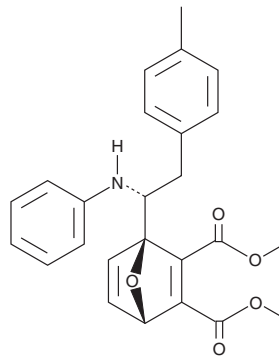
PRODUCT INFORMATION



COH000

Item No. 34906

CAS Registry No.: 1534358-79-6
Formal Name: (1R,4S)-1-[(1R)-2-(4-methylphenyl)-1-(phenylamino)ethyl]-7-oxabicyclo[2.2.1]hepta-2,5-diene-2,3-dicarboxylic acid, 2,3-dimethyl ester
MF: C₂₅H₂₅NO₅
FW: 419.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244 nm
Supplied as: A 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

COH000 is supplied as a solid. A stock solution may be made by dissolving the COH000 in the solvent of choice, which should be purged with an inert gas. COH000 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of COH000 in these solvents is approximately 13 and 11 mg/ml, respectively. It is also slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of COH000 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of COH000 in PBS (pH 7.2) is approximately 0.04 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

COH000 is an inhibitor of SUMO-activating enzyme (SAE; IC₅₀ = 0.2 μM), a heterodimer composed of SAE1 and SAE2.¹ It is selective for SAE over ubiquitin-activating enzyme (UAE) and NEDD8-activating enzyme (NAE) at 10 μM. COH000 induces degradation of SAE2 in HCT116 human colon cancer cells. COH000 (10 mg/kg) induces intratumoral apoptosis, decreases intratumoral c-Myc expression, and reduces tumor growth in an HCT116 mouse xenograft model.

Reference

1. Li, Y.-J., Du, L., Wang, J., et al. Allosteric inhibition of ubiquitin-like modifications by a class of inhibitor of SUMO-activating enzyme. *Cell Chem. Biol.* **26**(2), 278-288 (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, 11/15/2022

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