

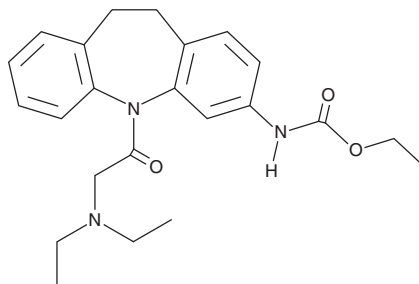
PRODUCT INFORMATION



CINPA1

Item No. 29110

CAS Registry No.: 102636-74-8
Formal Name: [5-[(diethylamino)acetyl]-10,11-dihydro-5H-dibenz[b,f]azepin-3-yl]-carbamic acid, ethyl ester
MF: C₂₃H₂₉N₃O₃
FW: 395.5
Purity: ≥98%
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

CINPA1 is supplied as a crystalline solid. A stock solution may be made by dissolving the CINPA1 in the solvent of choice, which should be purged with an inert gas. CINPA1 is soluble in organic solvents such as ethanol and DMSO. The solubility of CINPA1 in these solvents is approximately 100 mM.

Description

CINPA1 is a constitutive androstane receptor (CAR) antagonist (IC₅₀ = 70 nM in a reporter assay).¹ It is selective for CAR over the nuclear receptors GR, FXR, LXR α , LXR β , PPAR γ , RXR α , and RXR β at 18 μ M, but also antagonizes the human pregnane X receptor (PXR; IC₅₀ = 6.6 μ M). CINPA1 (0.3-5 μ M) inhibits CAR transactivation of CYP2B6 induced by CITCO (Item No. 16027) in primary human hepatocytes.

Reference

1. Cherian, M.T., Lin, W., Wu, J., *et al.* CINPA1 is an inhibitor of constitutive androstane receptor that does not activate pregnane X receptor. *Mol. Pharmacol.* **87**(5), 878-889 (2015).

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, 11/04/2022

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