

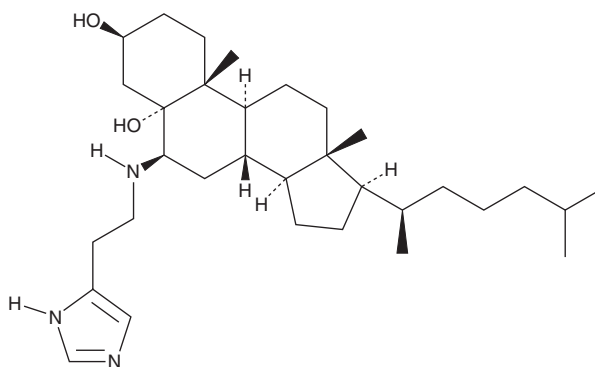
PRODUCT INFORMATION



Dendrogenin A

Item No. 37975

CAS Registry No.: 1191043-85-2
Formal Name: (3 β ,5 α ,6 β)-6-[[2-(1H-imidazol-5-yl)ethyl]amino]-cholestane-3,5-diol
Synonym: DDA
MF: C₃₂H₅₅N₃O₂
FW: 513.8
Purity: $\geq 95\%$
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

Dendrogenin A (DDA) is supplied as a solid. A stock solution may be made by dissolving the DDA in the solvent of choice, which should be purged with an inert gas. DDA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of DDA in these solvents is approximately 30 mg/ml.

Description

DDA is a selective liver X receptor (LXR) modulator (SLiM), an inhibitor of cholesterol epoxide hydrolase (ChEH; K_i = 120 nM), and an active metabolite of cholesterol.^{1,2} It is formed from 5,6 α -epoxy cholesterol via conjugation with histamine by DDA synthase.¹⁻³ DDA is found in non-cancerous human mammary epithelial cells and epithelial melanocytes but not in a variety of breast carcinoma or melanoma cells and only at low levels in isolated human breast tumor tissue.¹ It inhibits 22(R)-hydroxy cholesterol-induced activation of LXR β and LXR α in a reporter assay (IC₅₀s = 76 and 362 nM, respectively) but is also a partial agonist of LXRs, increasing protein levels of Nur77, NOR-1, LC3-I, and LC3-II in B16/F10 murine melanoma cells.⁴ It is selective for modulation of LXR α and LXR β over the pregnane X receptor (PXR), aryl hydrocarbon receptor (AhR), vitamin D receptor (VDR), retinoid X receptor γ (RXR γ), retinoic acid receptor α (RAR α), peroxisome proliferator-activated receptor α (PPAR α), PPAR γ , glucocorticoid receptor, androgen receptor, estrogen receptor α (ER α), and ER β at 2.5 μ M. It also increases protein levels of LC3-II in B16/F10 and SK-MEL-28 cancer cells when used at concentrations of 2.5 and 5 μ M and induces autophagic cell death in the same cell types at 2.5 μ M. DDA (0.37 μ g/kg) reduces tumor growth in a B16/F10 murine model of melanoma and a TS/A murine mammary cancer model and induces cancer cell differentiation *in vitro* and *in vivo*.²

References

- Poirot, M. and Silvente-Poirot, S. The tumor-suppressor cholesterol metabolite, dendrogenin A, is a new class of LXR modulator activating lethal autophagy in cancers. *Biochem. Pharmacol.* **153**, 75-81 (2018).
- de Medina, P., Paillasse, M.R., Segala, G., et al. Dendrogenin A arises from cholesterol and histamine metabolism and shows cell differentiation and anti-tumour properties. *Nat. Commun.* **4**, 1840 (2013).
- Soulès, R., Audouard-Combe, F., Huc-Claustre, E., et al. A fast UPLC-HILIC method for an accurate quantification of dendrogenin A in human tissues. *J. Steroid Biochem. Mol. Biol.* **194**, 105447 (2019).
- Segala, G., David, M., de Medina, P., et al. Dendrogenin A drives LXR to trigger lethal autophagy in cancers. *Nat. Commun.* **8**(1), 1903 (2017).

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, 03/27/2025

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