

PRODUCT INFORMATION



Vialinin A

Item No. 10010519

CAS Registry No.: 858134-23-3

Formal Name: benzeneacetic acid,
1,1'-(4,4'',5',6'-tetrahydroxy[1,1':4',1''-
terphenyl]-2',3'-diyl) ester

Synonym: Terrestrin A

MF: C₃₄H₂₆O₈

FW: 562.6

Purity: ≥95%

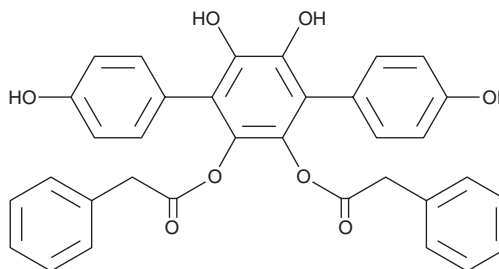
UV/Vis.: λ_{max}: 263 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Synthetic



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

Laboratory Procedures

Vialinin A is supplied crystalline solid. A stock solution may be made by dissolving the vialinin A in the solvent of choice, which should be purged with an inert gas. Vialinin A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of vialinin A in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Vialinin A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vialinin A should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Vialinin A has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Vialinin A is a terphenyl compound originally isolated from the fungi *T. terrestris* and *T. vialis*.^{1,2} Terphenyls, in general, are recognized as strong antioxidants.³ Vialinin A potently inhibits the release of TNF-α (IC₅₀ = 0.09 nM) and IL-4 (IC₅₀ = 2.8 nM), as well as β-hexosaminidase and CCL2 (MCP-1) from IgE-stimulated RBL-2H3 mast cells.⁴ Vialinin A does not significantly increase lactate dehydrogenase release from RBL-2H3 mast cells, suggesting low cytotoxicity.⁴

References

1. Radulovic, N., Quang, D.N., Hashimoto, T., et al. Terrestrins A-G: *p*-teraphenyl derivatives from the inedible mushroom *Thelephora terrestris*. *Phytochemistry* **66**(9), 1052-1059 (2005).
2. Xie, C., Koshino, H., Esumi, Y., et al. Vialinin A, a novel 2,2-diphenyl-1-picrylhydrazyl (DPPH) radical scavenger from an edible mushroom in China. *Biosci. Biotech. Biochem.* **69**(12), 2326-2332 (2005).
3. Liu, J.-K. Natural terphenyls: Developments since 1877. *Chem. Rev.* **106**(6), 2209-2223 (2006).
4. Onose, J., Xie, C., Ye, Y.Q., et al. Vialinin A, a novel potent inhibitor of TNF-α production from RBL-2H3 cells. *Biol. Pharm. Bull.* **31**(5), 831-833 (2008).

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, 01/29/2024

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