

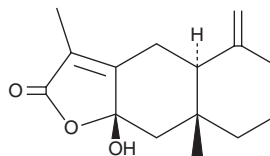
PRODUCT INFORMATION



Atractylenolide III

Item No. 24911

CAS Registry No.: 73030-71-4
Formal Name: (4aS,8aR,9aS)-4a,5,6,7,8,8a,9,9a-octahydro-9a-hydroxy-3,8a-dimethyl-5-methylene-naphtho[2,3-b]furan-2(4H)-one
Synonyms: Atractylenolide β , Codonolactone
MF: C₁₅H₂₀O₃
FW: 248.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 219 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Atractylenolide III is supplied as a crystalline solid. A stock solution may be made by dissolving the atractylenolide III in the solvent of choice. Atractylenolide III is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of atractylenolide III in ethanol is approximately 1 mg/ml and approximately 2 mg/ml in DMSO and DMF.

Atractylenolide III is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, atractylenolide III should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Atractylenolide III has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Atractylenolide III is a sesquiterpene that has been isolated from *Atractylodes* and has diverse biological activities, including anti-inflammatory, anti-angiogenic, pro- and anti-apoptotic, and neuroprotective properties.¹⁻⁵ It inhibits the release of nitric oxide (NO), prostaglandin E₂ (PGE₂; Item No. 14010), TNF- α , and IL-6 induced by LPS from RAW264.7 macrophages when used at a concentration of 50 μ M.¹ Atractylenolide III (100 μ M) induces apoptosis and cell cycle arrest, as well as increases caspase-3, caspase-9, and poly(ADP-ribose) polymerase (PARP) cleavage in A549 human lung carcinoma cells.³ However, in primary embryonic mouse cerebral cortical neurons, it inhibits glutamate-induced apoptosis and DNA fragmentation when used at a concentration of 40 μ M.⁴ It inhibits angiogenesis in an MDA-MB-231 human breast cancer mouse xenograft model when administered at a dose of 25 mg/kg per day.² Atractylenolide III (0.6 mg/kg per day) also decreases the latency to find the platform in the Morris water maze in a rat model of learning deficits induced by high-dose homocysteine.⁵

References

1. Ji, G.-Q., Chen, R.-Q., and Wang, L. *Immunopharmacol. Immunotoxicol.* **38**(2), 98-102 (2016).
2. Wang, S., Cai, R., Ma, J., et al. *Phytomedicine* **22**(11), 1017-1026 (2015).
3. Kang, T.-H., Bang, J.-Y., Kim, M.-H., et al. *Food Chem. Toxicol.* **49**(2), 514-519 (2011).
4. Liu, C., Zhao, H., Ji, Z.-H., et al. *Neurochem. Res.* **39**(9), 1753-1758 (2014).
5. Zhao, H., Ji, Z.-H., Liu, C., et al. *Neuroscience* **290**, 485-491 (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, 12/12/2022

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