

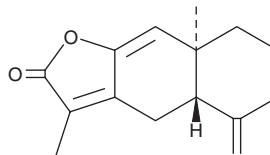
PRODUCT INFORMATION



Atractylenolide I

Item No. 25061

CAS Registry No.: 73069-13-3
Formal Name: 4aS,5,6,7,8,8aS-hexahydro-3,8a-dimethyl-5-methylene-naphtho[2,3-b]furan-2(4H)-one
MF: C₁₅H₁₈O₂
FW: 230.3
Purity: ≥98%
UV/Vis.: λ_{max}: 276 nm
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

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

Atractylenolide I is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, atractylenolide I should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Atractylenolide I 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 I is a sesquiterpene that has been found in the rhizomes of *A. macrocephala* and has diverse biological activities, including anti-inflammatory, anti-angiogenic, anti-tumor, and antidepressant properties.¹⁻⁴ It inhibits LPS-induced increases in TNF-α and nitric oxide (NO) production, as well as inducible nitric oxide synthase (iNOS) activity, in murine peritoneal macrophages (IC₅₀s = 23.1, 41, and 67.3 μM, respectively).¹ Atractylenolide I reduces pouch fluid weight, inflammatory cell count, granuloma weight, and vascular index (ID₅₀s = 24.18, 19.46, 14.44, and 15.15 mg/kg, respectively) in a mouse air pouch granuloma model induced by Freund's complete adjuvant (FCA).² It also reduces the number of microvessels in the air pouch wall by 58.27% when administered at a dose of 20 mg/kg and reduces the blood levels of TNF-α, IL-1β, IL-6, VEGF, placenta growth factor (PIGF), and bFGF in a dose-dependent manner in a mouse model of FCA-induced granuloma. Atractylenolide I inhibits the growth of T-24, 5637, RT4, and 253J bladder cancer cells *in vitro* (IC₅₀s = 12.8-63.7 μM) and reduces tumor growth in T-24 and 253J mouse xenograft models in a dose-dependent manner.³ In a mouse model of depression induced by chronic unpredictable mild stress, atractylenolide I reverses stress-induced decreases in hippocampal levels of serotonin (5-HT; Item No. 14332) and norepinephrine (Item No. 16673) and reduces immobility time in the forced swim and tail suspension tests in a dose-dependent manner, indicating antidepressant-like activity.⁴

References

1. Li, C.-Q., He, L.-C., and Jin, J.-Q. *Phytother. Res.* **21(4)**, 347-353 (2007).
2. Wang, C., Duan, H., and He, L. *Eur. J. Pharmacol.* **612(1-3)**, 143-152 (2009).
3. Yu, R., Yu, B.-X., Chen, J.-F., et al. *J. Exp. Clin. Cancer Res.* **35:40**, (2016).
4. Gao, H., Zhu, X., Xi, Y., et al. *Exp. Ther. Med.* **15(2)**, 1574-1579 (2018).

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, 10/21/2022

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