

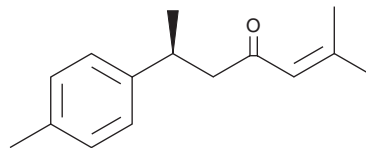
PRODUCT INFORMATION



(+)-ar-Turmerone

Item No. 11759

CAS Registry No.: 532-65-0
Formal Name: 2-methyl-6S-(4-methylphenyl)-2-hepten-4-one
Synonym: (S)-(+)-ar-Turmerone
MF: C₁₅H₂₀O
FW: 216.3
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

(+)-ar-Turmerone is an aromatic compound from the rhizomes of *C. longa*, a perennial plant of the ginger family, *Zingiberaceae*, commonly used in traditional Asian medicines.¹ It is also a major constituent of essential oil of turmeric, which is known for its antioxidant, anti-inflammatory, and anti-nociceptive properties.² Two structurally and physiologically distinct forms of turmerone have been isolated from *C. longa*, (+)-ar-Turmerone and α -turmerone.¹ (+)-ar-Turmerone stimulates the proliferation of peripheral blood mononuclear cells and increases the production of TNF- α , IL-2, and IFN- γ .¹ It also promotes the maturation of dendritic cells and induces the proliferation of neural stem cells both *in vitro* and *in vivo*.^{3,4}

References

1. Yue, G.G.L., Chan, B.C.L., Hon, P.M., *et al.* Evaluation of *in vitro* anti-proliferative and immunomodulatory activities of compounds isolated from *Curcuma longa*. *Food Chem. Toxicol.* **48(8-9)**, 2011-2020 (2010).
2. Liju, V.B., Jeena, K., and Kuttan, R. An evaluation of antioxidant, anti-inflammatory, and antinociceptive activities of essential oil from *Curcuma longa*. L. *Indian J. Pharmacol.* **43(5)**, 526-531 (2011).
3. Yonggang, T., Yiming, M., Heying, Z., *et al.* Maturation and upregulation of functions of murine dendritic cells (DCs) under the influence of purified aromatic-turmerone (AR). *Hum. Vaccin. Immunother.* **8(10)**, 1416-1424 (2012).
4. Hucklenbrioch, J., Klein, R., Neumaier, B., *et al.* Aromatic-turmerone induces neural stem cell proliferation *in vitro* and *in vivo*. *Stem Cell Res. Ther.* **5(4)**, 100 (2014).

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, 02/21/2023

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