

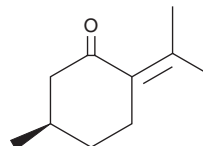
PRODUCT INFORMATION



(R)-(+)-Pulegone

Item No. 23163

CAS Registry No.: 89-82-7
Formal Name: 5R-methyl-2-(1-methylethylidene)-cyclohexanone
Synonyms: NSC 15334, D-Pulegone
MF: C₁₀H₁₆O
FW: 152.2
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(R)-(+)-Pulegone is supplied as a neat oil. A stock solution may be made by dissolving the (R)-(+)-pulegone in the solvent of choice. (R)-(+)-Pulegone is soluble in organic solvents such as ethanol and chloroform.

Description

(R)-(+)-Pulegone is a monoterpene that has been found in a variety of plants, including *Labiatae* species and *M. verticillata*, and has anti-inflammatory activity.¹ It stimulates proliferation of lymphocytes derived from allergic patients *in vitro* when used at concentrations of 6.2 and 62 µg/ml and reduces the level of IL-13 in supernatant from lymphocyte cultures when used at a concentration of 62 µg/ml.² (R)-(+)-Pulegone decreases scratching behavior and reduces skin thickness in a mouse model of atopic dermatitis when administered topically at doses ranging from 0.1 to 10 µM. It also decreases serum IgE levels and skin protein levels of nerve growth factor, IFN-γ, IL-6, IL-1β, TNF-α, NF-κB, phosphorylated ERK1/2, and phosphorylated JNK.

References

1. Choi, Y.Y., Kim, M.H., Lee, H., *et al.* (R)-(+)-pulegone suppresses allergic and inflammation responses on 2,4-dinitrochlorobenzene-induced atopic dermatitis in mice model. *J. Dermatol. Sci.* (2018).
2. Cariddi, L., Escobar, F., Moser, M., *et al.* Monoterpenes isolated from *Minthostachys verticillata* (Griseb.) Epling essential oil modulates immediate-type hypersensitivity responses *in vitro* and *in vivo*. *Planta Med.* **77**(15), 1687-1694 (2011).

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, 11/28/2022

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