

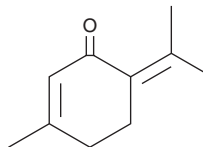
PRODUCT INFORMATION



Piperitenone

Item No. 25752

CAS Registry No.: 491-09-8
Formal Name: 3-methyl-6-(1-methylethylidene)-2-cyclohexen-1-one
Synonym: NSC 667470
MF: $C_{10}H_{14}O$
FW: 150.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 244, 278 nm
Supplied as: A 10 mg/ml solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Piperitenone is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Piperitenone is miscible in ethanol, DMSO, and dimethyl formamide.

Description

Piperitenone is a monoterpene that has been found in various plants, including *Cannabis*, and has antioxidant activity.^{1,2} It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in a cell free assay ($IC_{50} = 22.7 \mu g/ml$).² Piperitenone also inhibits peroxidation of linoleic acid (Item No. 90150).

References

1. Turner, C.E., Elsohly, M.A., and Boeren, E.G. Constituents of *Cannabis sativa* L. XVII. A review of the natural constituents. *J. Nat. Prod.* **43**(2), 169-234 (1980).
2. Iqbal, T., Hussain, A.I., Chatha, S.A., *et al.* Antioxidant activity and volatile and phenolic profiles of essential oil and different extracts of wild mint (*Mentha longifolia*) from the Pakistani flora. *J. Anal. Methods Chem.* **536490** (2013).

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, 03/25/2026

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