

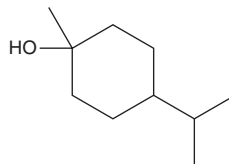
PRODUCT INFORMATION



4-Isopropyl-1-methylcyclohexanol

Item No. 45695

CAS Registry No.: 21129-27-1
Formal Name: 1-methyl-4-(1-methylethyl)-cyclohexanol
Synonym: *p*-Methan-1-ol
MF: C₁₀H₂₀O
FW: 456.3
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

4-Isopropyl-1-methylcyclohexanol is supplied as a liquid. A stock solution may be made by dissolving the 4-isopropyl-1-methylcyclohexanol in the solvent of choice, which should be purged with an inert gas. 4-Isopropyl-1-methylcyclohexanol is sparingly soluble (1-10 mg/ml) in ethanol.

Description

4-Isopropyl-1-methylcyclohexanol is a dihydroterpineol that has been found in *Mentha x piperita*.^{1,2}

References

1. Xu, S., Zhao, Z., Bi, L., *et al.* Selective aerobic hydroxylation of *p*-menthane to dihydroterpineols catalyzed by metalloporphyrins in solvent and additive free system. *Catal. Commun.* **59**, 26-29 (2015).
2. Yurchiak, M.E., Bailey, S., Sakib, A.H., *et al.* Aqueous leaf extracts of peppermint (*Mentha x piperita*) and white snakeroot (*Ageratina altissima*) exhibit antibacterial and antiviral activity. *Microorganisms* **14**(1), 80 (2025).

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, 06/11/2026

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