

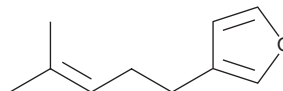
PRODUCT INFORMATION



Perillene

Item No. 35006

CAS Registry No.: 539-52-6
Formal Name: 3-(4-methyl-3-penten-1-yl)-furan
MF: C₁₀H₁₄O
FW: 150.2
Purity: ≥98%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Perilla frutescens*



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

Laboratory Procedures

Perillene is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of perillene in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Perillene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the acetonitrile solution of perillene should be diluted with the aqueous buffer of choice. Perillene has a solubility of 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Perillene is a monoterpene that has been found in *P. citriodora*.^{1,2}

References

1. Umney, J.C. Essential oil of perilla. *Perfum. Essent. Oil Rec.* **10**, 170-171 (1919).
2. Nishizawa, A., Honda, G., and Tabata, M. Genetic control of perillene accumulation in *Perilla frutescens*. *Phytochemistry* **29(9)**, 2873-2875 (1990).

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, 02/17/2026

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