

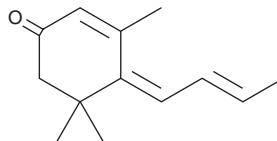
PRODUCT INFORMATION



Megastigmatrienone (mixed isomers)

Item No. 38754

CAS Registry No.: 13215-88-8
Formal Name: 4-(2-buten-1-ylidene)-3,5,5-trimethyl-2-cyclohexen-1-one
Synonym: Tabanone
MF: C₁₃H₁₈O
FW: 190.3
Supplied as: A neat 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

Megastigmatrienone (mixed isomers) is supplied as a neat liquid. A stock solution may be made by dissolving the megastigmatrienone (mixed isomers) in the solvent of choice, which should be purged with an inert gas. Megastigmatrienone (mixed isomers) is soluble (≥10 mg/ml) in ethanol and DMSO and slightly soluble (0.1-1 mg/ml) in methanol.

Description

Megastigmatrienone (mixed isomers) is a mixture of isomers of megastigmatrienone, a compound that has the aroma of tobacco.¹ Isomers of megastigmatrienone have been found in tobacco and oak barrel-aged wine.^{1,2}

References

1. Fujimori, T., Kasuga, R., Matsushita, H., *et al.* Neutral volatile components of burley tobacco. *Agr. Biol. Chem.* **40**(2), 303-315 (1976).
2. Slaghenaufl, D., Perello, M.-C., Marchand-Marion, S., *et al.* Quantitative solid phase microextraction--gas chromatography mass spectrometry analysis of five megastigmatrienone isomers in aged wine. *Anal. Chim. Acta* **813**, 63-69 (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, 07/23/2024

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