

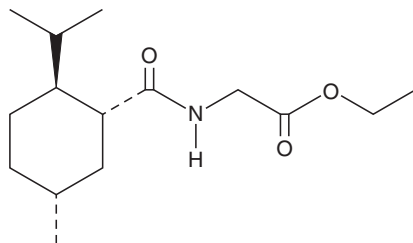
PRODUCT INFORMATION



WS5

Item No. 38763

CAS Registry No.: 68489-14-5
Formal Name: N-[[[(1R,2S,5R)-5-methyl-2-(1-methylethyl)cyclohexyl]carbonyl]-glycine, ethyl ester
MF: $C_{15}H_{27}NO_3$
FW: 269.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

WS5 is supplied as a solid. A stock solution may be made by dissolving the WS5 in the solvent of choice, which should be purged with an inert gas. WS5 is slightly soluble in chloroform and methanol.

Description

WS5 is a cooling compound and a derivative of the monoterpene (-)-menthol (Item No. 30384).¹ Formulations containing WS5 have been used as cooling agents in toothpaste and hard candies, as well as skin creams and lotions.

Reference

1. Watson, H.R., Hems, R., Rowsell, D.G., *et al.* New compounds with the menthol cooling effect. *J. Soc. Cosmet. Chem.* **29**, 185-200 (1978).

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, 12/07/2023

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