

# PRODUCT INFORMATION



## 2-Methylhexacosane

Item No. 9003585

CAS Registry No.: 1561-02-0

Synonym: 2-Me-C26

MF:  $C_{27}H_{56}$

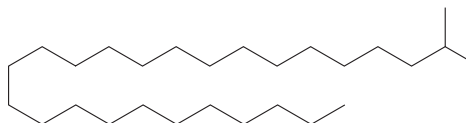
FW: 380.7

Purity:  $\geq 95\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

2-Methylhexacosane is supplied as a solid. A stock solution may be made by dissolving the 2-methylhexacosane in the solvent of choice, which should be purged with an inert gas. 2-Methylhexacosane is soluble in chloroform and hexane.

### Description

2-Methylhexacosane is a saturated hydrocarbon and an insect pheromone.<sup>1,2</sup> It has been found in the cuticle of *M. dasystomus* females, but not males, where it contributes to the mating behavior of males, as well as in *D. melanogaster* females where it modulates aggression of males towards females. 2-Methylhexacosane has also been found in yellow jacket (*V. vulgaris*) trail extracts.<sup>3</sup>

### References

1. Spikes, A.E., Pashen, M.A., Millar, J.G., *et al.* First contact pheromone identified for a longhorned beetle (Coleoptera: Cerambycidae) in the subfamily Prioninae. *J. Chem. Ecol.* **36**(9), 943-954 (2010).
2. Fernández, d.I.P., Chan, Y.-B., Yew, J.Y., *et al.* Pheromonal and behavioral cues trigger male-to-female aggression in *Drosophila*. *PLoS Biol.* **8**(11), e1000541 (2010).
3. Steinmetz, I., Schmolz, E., and Ruther, J. Cuticular lipids as trail pheromone in a social wasp. *Proc. Biol. Sci.* **270**(1513), 385-391 (2003).

#### 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, 12/16/2022

#### 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