

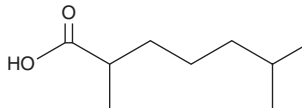
PRODUCT INFORMATION



DL-2,6-dimethyl Heptanoic Acid

Item No. 24822

CAS Registry No.: 7494-12-4
Formal Name: 2,6-dimethyl-heptanoic acid
Synonyms: 2,6-Dimethyl C7:0 Fatty Acid,
FA 9:0
MF: C₉H₁₈O₂
FW: 158.2
Purity: ≥98%
Supplied as: A solution in chloroform
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

DL-2,6-dimethyl Heptanoic acid is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. DL-2,6-dimethyl Heptanoic acid is soluble in the organic solvent chloroform.

Description

DL-2,6-dimethyl Heptanoic acid is an aliphatic branched acid that inhibits the growth of *V. fischeri* (EC₅₀ = 17 μM).¹ It is a building block in the synthesis of 2,6-dimethyl heptanoyl-CoA, a substrate of acyl-CoA dehydrogenase.²

References

1. Jones, D., Scarlett, A.G., West, C.E., *et al.* Toxicity of individual naphthenic acids to *Vibrio fischeri*. *Environ. Sci. Technol.* **45**(22), 9776-9782 (2011).
2. Sobhi, H.F., Minkler, P.E., and Hoppel, C.L. Synthesis and characterization of *cis*-4-decenoyl-CoA, 3-phenylpropionyl-CoA, and 2,6-dimethylheptanoyl-CoA. *Anal. Biochem.* **401**(1), 114-124 (2010).

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, 03/06/2024

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