

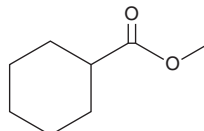
PRODUCT INFORMATION



Methyl cyclohexanecarboxylate

Item No. 30403

CAS Registry No.: 4630-82-4
Formal Name: cyclohexanecarboxylic acid, methyl ester
MF: $C_8H_{14}O_2$
FW: 142.2
Purity: $\geq 95\%$
Supplied as: A liquid
Storage: $-20^\circ C$
Stability: ≥ 2 years



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

Laboratory Procedures

Methyl cyclohexanecarboxylate is supplied as a liquid. A stock solution may be made by dissolving the methyl cyclohexanecarboxylate in the solvent of choice, which should be purged with an inert gas. Methyl cyclohexanecarboxylate is soluble in ethanol, DMSO, and dimethyl formamide. The solubility of methyl cyclohexanecarboxylate in these solvents is approximately 30 mg/ml.

Description

Methyl cyclohexanecarboxylate is an aliphatic ester that has been used in a variety of chemical synthesis studies, including the preparation of primary amides and reduction of esters to alcohols.^{1,2}

References

1. Prasanth, C.P., Ebbin, J., Abhijith, A., *et al.* Stabilization of $NaBH_4$ in methanol using a catalytic amount of NaOMe. Reduction of esters and lactones at room temperature without solvent-induced loss of hydride. *J. Org. Chem.* **83**(3), 1431-1440 (2018).
2. Veitch, G.E., Bridgwood, K.L., and Ley, S.V. Magnesium nitride as a convenient source of ammonia: Preparation of primary amides. *Org. Lett.* **10**(16), 3623-3625 (2008).

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, 05/11/2020

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