

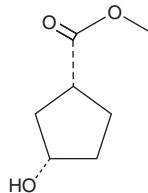
PRODUCT INFORMATION



(1R,3S)-3-Hydroxycyclopentane carboxylic acid methyl ester

Item No. 11374

CAS Registry No.: 174292-59-2
Formal Name: (1R,3S)-3-hydroxy-cyclopentanecarboxylic acid, methyl ester
MF: C₇H₁₂O₃
FW: 144.2
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(1R,3S)-3-Hydroxycyclopentane carboxylic acid methyl ester is supplied as a neat oil. A stock solution may be made by dissolving the (1R,3S)-3-hydroxycyclopentane carboxylic acid methyl ester in the solvent of choice, which should be purged with an inert gas. (1R,3S)-3-Hydroxycyclopentane carboxylic acid methyl ester is soluble in the organic solvent DMSO.

Description

(1R,3S)-3-Hydroxycyclopentane carboxylic acid methyl ester is a synthetic intermediate useful for pharmaceutical synthesis.

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, 12/12/2022

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