

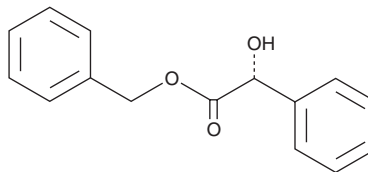
PRODUCT INFORMATION



(R)-benzyl Mandelate

Catalog No. 10007145

CAS Registry No.: 97415-09-3
Formal Name: α-hydroxy-benzeneacetic acid, phenylmethyl ester
Synonyms: Benzyl D-mandelate
MF: C₁₅H₁₄O₃
FW: 242.3
Purity: ≥99%
Supplied as: A crystalline 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

(R)-benzyl Mandelate is supplied as a crystalline solid. A stock solution may be made by dissolving the (R)-benzyl mandelate in the solvent of choice, which should be purged with an inert gas. (R)-benzyl Mandelate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (R)-benzyl mandelate in these solvents is approximately 20, 25, and 30 mg/ml, respectively.

(R)-benzyl Mandelate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (R)-benzyl mandelate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (R)-benzyl Mandelate has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(R)-benzyl Mandelate 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, 11/16/2022

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