

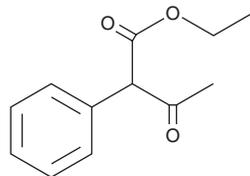
PRODUCT INFORMATION



Ethyl 2-phenylacetoacetate

Item No. 39487

CAS Registry No.: 5413-05-8
Formal Name: α -acetyl-benzeneacetic acid, ethyl ester
Synonyms: EAPA, Ethyl α -acetylbenzeneacetate, Ethyl α -acetylphenylacetate, Ethyl α -phenylacetoacetate, Ethyl 3-oxo-2-phenylbutanoate, Ethyl 3-oxo-2-phenylbutyrate, NSC 6389



MF: $C_{12}H_{14}O_3$
FW: 206.2
Purity: $\geq 98\%$ (mixture of isomers)
Supplied as: A neat oil
Storage: $-20^{\circ}C$
Stability: ≥ 2 years

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

Description

Ethyl 2-phenylacetoacetate (Item No. 39487) is an analytical reference standard categorized as a precursor in the synthesis of the amphetamine precursor phenylacetone (Item Nos. 16103 | 16444).¹ This product is intended for research and forensic applications.

Reference

1. Tsujikawa, K., Okada, Y., Segawa, H., *et al.* Analysis of potential phenylacetone precursors (ethyl 3-oxo-2-phenylbutyrate, methyl 3-oxo-4-phenylbutyrate, and ethyl 3-oxo-4-phenylbutyrate) by gas chromatography/mass spectrometry and their conversion to phenylacetone. *Drug Test Anal.* **14**(3), 439-449 (2022).

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, 02/19/2024

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