

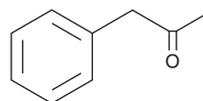
PRODUCT INFORMATION



Phenylacetone (exempt preparation)

Item No. 16444

CAS Registry No.: 103-79-7
Formal Name: 1-phenyl-2-propanone
Synonyms: Benzyl Methyl Ketone, P2P, Phenyl-2-propanone
MF: C₉H₁₀O
FW: 134.2
Purity: ≥98%
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥3 years



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

Description

Phenylacetone (exempt preparation) (Item No. 16444) is an analytical reference standard categorized as a precursor in the synthesis of amphetamine and methamphetamine.^{1,2} Phenylacetone is regulated as a Schedule II compound in the United States. Phenylacetone (exempt preparation) (Item No. 16444) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Krawczyk, W., Kunda, T., Perkowska, I., *et al.* Impurity profiling/comparative analyses of samples of 1-phenyl-2-propanone. *Bull. Narc.* **57(1-2)**, 33-62 (2005).
2. Sinnema, A. and Verweij, A.M.A. Impurities in illicit amphetamine: A review. *Bull. Narc.* **33(3)**, 37-54 (1981).

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, 01/15/2025

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