

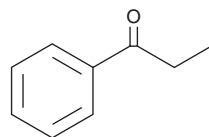
# PRODUCT INFORMATION



## Propiophenone

Item No. 41904

CAS Registry No.: 93-55-0  
Formal Name: 1-phenyl-1-propanone  
Synonyms: NSC 16937, Phenyl ethyl ketone, Propionylbenzene  
MF:  $C_9H_{10}O$   
FW: 134.2  
Purity:  $\geq 98\%$   
Supplied as: A 50 mg/ml solution in acetonitrile  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 1$  year



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

### Description

Propiophenone (Item No. 41904) is an analytical reference standard categorized as a precursor in the synthesis of cathinones.<sup>1</sup> This product is intended for research and forensic applications.

### Reference

1. Huxley, C.M., Biddle, T., Shand, E., *et al.* Forensic implications of novel synthesis of cathinone derivatives by Neber and modified Neber rearrangements. *Forensic Chem.* **38(6602)**, 100558 (2024).

#### 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, 09/18/2025

#### 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