

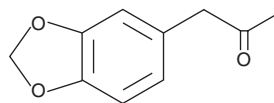
PRODUCT INFORMATION



Piperonyl methyl ketone

Item No. 13892

CAS Registry No.: 4676-39-5
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-propanone
Synonyms: MDP2P, 3,4-Methylenedioxyphenylpropan-2-one, NSC 16688, PMK, PVI-2
MF: C₁₀H₁₀O₃
FW: 178.2
Purity: ≥95%
UV/Vis.: λ_{max}: 237, 287 nm
Supplied as: A 50 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

Piperonyl methyl ketone (Item No. 13892) is an analytical reference standard categorized as a precursor in the synthesis of methylenedioxy phenethylamines and amphetamines, including 3,4-MDMA (Item Nos. ISO60190 | 13971).¹ Piperonyl methyl ketone is also a potential impurity found in illicit preparations of methylenedioxy phenethylamines.² This product is intended for research and forensic applications.

References

- Schäffer, M., Dieckmann, S., Pütz, M., *et al.* Impact of reaction parameters on the chemical profile of 3,4-methylenedioxymethamphetamine synthesized *via* reductive amination: Target analysis based on GC-qMS compared to non-targeted analysis based on GC × GC-TOF-MS. *Forensic Sci. Int.* **233(1-3)**, 201-211 (2013).
- Swist, M., Wilamowski, J., and Parczewski, A. Determination of synthesis method of ecstasy based on the basic impurities. *Forensic Sci. Int.* **152(2-3)**, 175-184 (2005).

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, 12/05/2025

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