

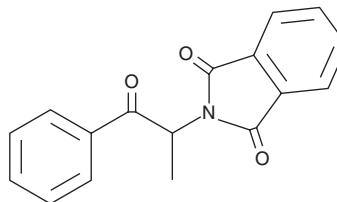
PRODUCT INFORMATION



α -Phthalimidopropiophenone

Item No. 17141

CAS Registry No.: 19437-20-8
Formal Name: 2-(1-methyl-2-oxo-2-phenylethyl)-1H-isoindole-1,3(2H)-dione
Synonyms: α -PAPP,
2-Phthalimidopropiophenone
MF: C₁₇H₁₃NO₃
FW: 279.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 208, 221, 291 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

α -Phthalimidopropiophenone is a cathinone analog that has been identified in capsules being distributed by the illicit drug market.¹ The substitution of the phthalimido moiety for the typical α -amino group may increase the stability of this cathinone in storage. Alternatively, similar nitrogen-protected propiophenones have been shown to be metabolized *in vivo* to produce cathinone, suggesting that this compound may act as a prodrug.^{1,2} The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

References

1. Camilleri, A., Johnston, M.R., Brennan, M., *et al.* Chemical analysis of four capsules containing the controlled substance analogues 4-methylmethcathinone, 2-fluoromethamphetamine, α -phthalimidopropiophenone and N-ethylcathinone. *Forensic Sci. Int.* **197(1-3)**, 59-66 (2010).
2. Springer, D., Fritschi, G., and Maurer, H. Metabolism of the new designer drug α -pyrrolidinopropiophenone (PPP) and the toxicological detection of PPP and 4'-methyl- α -pyrrolidinopropiophenone (MPPP) studied in rat urine using gas chromatography-mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **796(2)**, 253-266 (2003).

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, 07/05/2023

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