

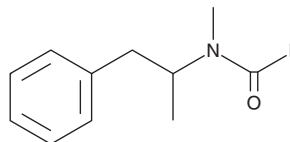
PRODUCT INFORMATION



N-Formylmethamphetamine

Item No. 15634

CAS Registry No.: 42932-20-7
Formal Name: N-methyl-N-(1-methyl-2-phenylethyl)-formamide
MF: $C_{11}H_{15}NO$
FW: 177.2
Purity: $\geq 98\%$
Supplied as: A solution in methyl acetate
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Description

N-Formylmethamphetamine is an impurity generated in the synthesis of methamphetamine.^{1,2} While suggested as being a route-specific impurity of the Leuckart method, this impurity has also been detected in methamphetamine samples produced from ephedrine or by reductive amination.^{3,4} The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

References

1. Kuwayama, K., Tsujikawa, K., Miyaguchi, H., *et al.* Identification of impurities and the statistical classification of methamphetamine using headspace solid phase microextraction and gas chromatography-mass spectrometry. *Forensic Sci. Int.* **160**, 44-52 (2006).
2. Qi, Y., Evans, I.D., and McCluskey, A. Australian Federal Police seizures of illicit crystalline methamphetamine ('ice') 1998-2002: Impurity analysis. *Forensic Sci. Int.* **164**, 201-210 (2006).
3. Qi, Y., Evans, I., and McCluskey, A. New impurity profiles of recent Australian imported 'ice': Methamphetamine impurity profiling and the identification of (pseudo)ephedrine and Leuckart specific marker compounds. *Forensic Sci. Int.* **169(2-3)**, 173-180 (2007).
4. Kunalan, V., Nic Daéid, N., Kerr, W.J., *et al.* Characterization of route specific impurities found in methamphetamine synthesized by the Leuckart and reductive amination methods. *Anal. Chem.* **81(17)**, 7342-7348 (2009).

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, 06/23/2023

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