

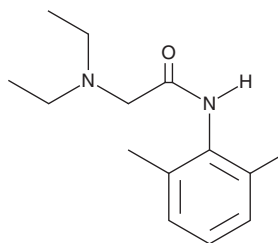
PRODUCT INFORMATION



Lidocaine

Item No. 20081

CAS Registry No.: 137-58-6
Formal Name: 2-(diethylamino)-N-(2,6-dimethylphenyl)-acetamide
Synonym: NSC 40030
MF: C₁₄H₂₂N₂O
FW: 234.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Lidocaine (Item No. 20081) is an analytical reference standard that is categorized as an adulterant. It is commonly used as a topical anesthetic.¹ Lidocaine is also found as an adulterant in cocaine.²⁻⁵ This product is intended for research and forensic applications.

References

1. Kumar, M., Chawla, R., and Goyal, M., Topical anesthesia. *J. Anaesthesiol. Clin. Pharmacol.* **31(4)**, 450-456 (2015).
2. DeCarvalho, T.C., Tosato, F., Souza, L.M., *et al.* Thin layer chromatography coupled to paper spray ionization mass spectrometry for cocaine and its adulterants analysis. *Forensic Sci. Int.* **262**, 56-65 (2016).
3. Floriani, G., Gasparetto, J.C., Pontarolo, R., *et al.* Development and validation of an HPLC-DAD method for simultaneous determination of cocaine, benzoic acid, benzoylecgonine and the main adulterants found in products based on cocaine. *Forensic Sci. Int.* **235**, 32-39 (2014).
4. Grobério, T.S., Zacca, J.J., Botelho, É.D., *et al.* Discrimination and quantification of cocaine and adulterants in seized drug samples by infrared spectroscopy and PLSR. *Forensic Sci. Int.* **257**, 297-306 (2015).
5. Phillips, K.A., Hirsch, G.A., Epstein, D.H., *et al.* Cardiac complications of unwitting co-injection of quinine/quinidine with heroin in an intravenous drug user. *J. Gen. Intern. Med.* **27(12)**, 1722-1725 (2012).

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, 02/29/2024

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