

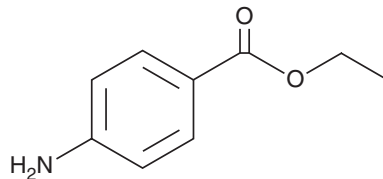
# PRODUCT INFORMATION



## Benzocaine

Item No. 20132

**CAS Registry No.:** 94-09-7  
**Formal Name:** 4-amino-benzoic acid, ethyl ester  
**Synonyms:** Ethyl 4-Aminobenzoate, H-4-Abz-OEt, NSC 4688, NSC 41531  
**MF:** C<sub>9</sub>H<sub>11</sub>NO<sub>2</sub>  
**FW:** 165.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 294 nm  
**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

Benzocaine (Item No. 20132) is an analytical reference standard that is categorized as an adulterant. It is commonly used as a topical anesthetic.<sup>1</sup> Benzocaine is also found as an adulterant in cocaine.<sup>2-5</sup> 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., L. 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