

PRODUCT INFORMATION



Cocaine (hydrochloride)

Item No. 33764

CAS Registry No.: 53-21-4
Formal Name: (1R,2R,3S,5S)-3-(benzoyloxy)-8-(methyl)-8-azabicyclo[3.2.1]octane-2-carboxylic acid, methyl ester, monohydrochloride

Synonyms: Benzoylmethylecgonine, L-Cocaine, NSC 25263

MF: C₁₇H₂₁NO₄ • HCl

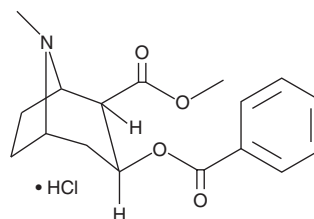
FW: 339.8

Purity: ≥98%

Supplied as: A neat solid

Storage: -20°C

Stability: ≥7 years



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

Description

Cocaine (hydrochloride) (Item No. 33764) is an analytical reference material categorized as a tropane. Cocaine has a high potential for abuse and has been found in samples seized by law enforcement.^{1,2} Cocaine is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Hemby, S.E. Cocainomics: New insights into the molecular basis of cocaine addiction. *J. Neuroimmune Pharmacol.* **5**(1), 70-82 (2010).
2. 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).

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, 06/22/2021

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