

PRODUCT INFORMATION



Norcocaethylene (hydrochloride)

Item No. 39838

CAS Registry No.: 362044-01-7

Formal Name: (1R,2R,3S,5S)-3-(benzoyloxy)-8-azabicyclo[3.2.1]octane-2-carboxylic acid, monohydrochloride, ethyl ester

Synonyms: Benzoylnorecgonine ethyl ester, N-Demethylcocaethylene, Norethylcocaine, Norhomococaine

MF: $C_{17}H_{21}NO_4 \cdot HCl$

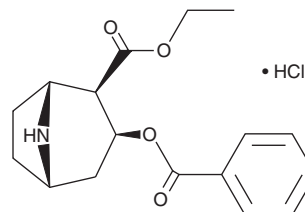
FW: 339.8

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Description

Norcocaethylene (hydrochloride) (Item No. 39838) is an analytical reference standard categorized as a tropane metabolite.¹ Norcocaethylene is a toxic metabolite of cocaine (Item Nos. ISO60176 | 16186 | 33764) and cocaethylene (Item No. 28820) that occurs with concurrent use of cocaine and alcohol.^{1,2} Norcocaethylene is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

References

1. Dean, R.A., Harper, E.T., Dumaul, N., *et al.* Effects of ethanol on cocaine metabolism: Formation of cocaethylene and norcocaethylene. *Toxicol. Appl. Pharmacol.* **117**(1), 1-8 (1992).
2. Zheng, X., Shang, L., Zhan, C.G., *et al.* In vivo characterization of toxicity of norcocaethylene and norcocaine identified as the most toxic cocaine metabolites in male mice. *Drug Alcohol Depend.* **204**, 107462 (2019).

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, 02/11/2025

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