

PRODUCT INFORMATION



Cocaethylene-d₃

Item No. 18038

CAS Registry No.: 136765-30-5

Formal Name: (1R,2R,3S,5S)-3-(benzoyloxy)-8-(methyl-d₃)-8-azabicyclo[3.2.1]octane-2-carboxylic acid, ethyl ester

Synonyms: Ethyl benzoylecgonine-d₃, Ethylcocaine-d₃, Homocaine-d₃

MF: C₁₈H₂₀D₃NO₄

FW: 320.4

Chemical Purity: ≥98% (cocaethylene)

Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

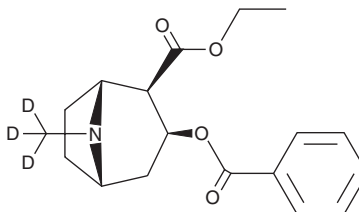
Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Special conditions: No protic solvents - will transesterify

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



Description

Cocaethylene-d₃ (Item No. 18038) is intended for use as an internal standard for the quantification of cocaethylene (Item No. 28820) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Cocaethylene is categorized as a metabolite of cocaine in the presence of ethanol.¹ It has been found in the blood of individuals following ingestion of cocaine concomitantly with ethanol, and it has reinforcing properties in rhesus monkeys. This product is intended for research and forensic applications.

Reference

1. Jatlow, P., Elsworth, J.D., Bradberry, C.W., et al. Cocaethylene: A neuropharmacologically active metabolite associated with concurrent cocaine-ethanol ingestion. *Life Sci.* **48(18)**, 1787-1794 (1991).

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, 01/27/2025

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