

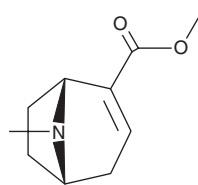
PRODUCT INFORMATION



Anhydroecgonine methyl ester

Item No. 17979

CAS Registry No.: 43021-26-7
Formal Name: (1R,5S)-8-methyl-8-azabicyclo[3.2.1]oct-2-ene-2-carboxylic acid, methyl ester
Synonyms: AEME, Methylecgonidine
MF: $C_{10}H_{15}NO_2$
FW: 181.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 218 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

Anhydroecgonine methyl ester (Item No. 17979) is an analytical reference standard categorized as a pyrolysis product of cocaine.¹ Anhydroecgonine methyl ester has been used as a biomarker for smoked cocaine. Anhydroecgonine methyl ester is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

Reference

1. Riley, K.J., Lu, N.T., Meeker, J.E., *et al.* Monitoring the crack epidemic through urine testing: Establishment of routine detection methods. *Addict. Biol.* **6**, 83-95 (2001).

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, 09/18/2025

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