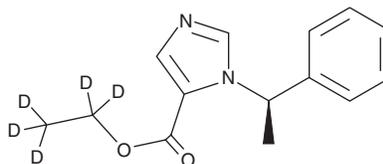


PRODUCT INFORMATION



Etomidate-d₅ Item No. 44952

CAS Registry No.: 3046414-07-4
Formal Name: 1-[(1R)-1-phenylethyl]-1H-imidazole-5-carboxylic acid ethyl-d₅ ester
Synonyms: (+)-Etomidate-d₅, d-Etomidate-d₅, (R)-Etomidate-d₅
MF: C₁₄H₁₁D₅N₂O₂
FW: 249.3
Chemical Purity: ≥98% (Etomidate)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

Etomidate-d₅ (Item No. 44952) is an analytical reference standard intended for use as an internal standard for the quantification of etomidate (Item Nos. 41826 | 21993) 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).

Etomidate is categorized as an anesthetic.^{1,2} Etomidate has been associated with overdose and death.³ Formulations containing etomidate have been used as anesthetics. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 40297).

References

1. Brede, M., Philipp, M., Knaus, A., *et al.* α_2 -adrenergic receptor subtypes - novel functions uncovered in gene-targeted mouse models. *Biol. Cell* **96**(5), 343-348 (2004).
2. Green, C.J., Knight, J., Precious, S., *et al.* Metomidate, etomidate and fentanyl as injectable anaesthetic agents in mice. *Lab. Anim.* **15**(2), 171-175 (1981).
3. Molina, D.K., Hargrove, V.M., and Rodriguez, R.G. Distribution of etomidate in a fatal intoxication. *J. Anal. Toxicol.* **32**(8), 715-718 (2008).

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, 12/17/2025

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