

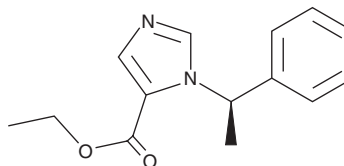
PRODUCT INFORMATION



Etomidate

Item No. 21993

CAS Registry No.: 33125-97-2
Formal Name: 1-[(1R)-1-phenylethyl]-1H-imidazole-5-carboxylic acid ethyl ester
Synonyms: (+)-Etomidate, d-Etomidate, (R)-Etomidate, R 16659
MF: C₁₄H₁₆N₂O₂
FW: 244.3
Purity: ≥98%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Etomidate (Item No. 21993) is an analytical reference standard that is categorized as an anesthetic. Formulations containing etomidate are used for general anesthesia and sedation.¹ Etomidate also inhibits adrenal cortisol biosynthesis and mitigates hypercortisolism, as is found in Cushing syndrome.^{2,3} This product is intended for research and forensic applications.

References

1. Forman, S.A. Clinical and molecular pharmacology of etomidate. *Anesthesiology* **114**(3), 695-707 (2011).
2. Juszczak, A. and Grossman, A. The management of Cushing's disease - from investigation to treatment. *Endokrynol. Pol.* **64**(2), 166-174 (2013).
3. Preda, V.A., Sen, J., Karavitaki, N., et al. Etomidate in the management of hypercortisolaemia in Cushing's syndrome: A review. *Eur. J. Endocrinol.* **167**(2), 137-143 (2012).

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, 04/01/2024

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