

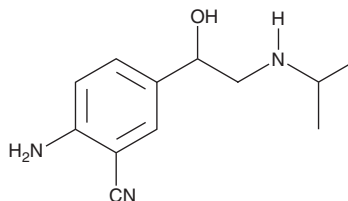
PRODUCT INFORMATION



Cimaterol

Item No. 15613

CAS Registry No.: 54239-37-1
Formal Name: 2-amino-5-[1-hydroxy-2-[(1-methylethyl)amino]ethyl]-benzonitrile
Synonym: Cimaterolum
MF: C₁₂H₁₇N₃O
FW: 219.3
Purity: ≥99%
UV/Vis.: λ_{max}: 218, 253 nm
Supplied as: A solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

Cimaterol is supplied as a solid. A stock solution may be made by dissolving the cimaterol in the solvent of choice. Cimaterol is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of cimaterol in these solvents is approximately 50 and 100 mM, respectively.

Description

Cimaterol is a potent agonist of β-adrenergic receptors (pEC₅₀ = 8.13, 8.78, and 6.62 for human β₁, β₂, and β₃, respectively).¹ Formulations containing cimaterol have been used in farmed animals (swine, fowl, etc.) to increase carcass mass and to alter muscle and fat deposition.^{2,3}

References

1. Mistry, S.N., Baker, J.G., Fischer, P.M., *et al.* Synthesis and in vitro and in vivo characterization of highly β₁-selective β-adrenoceptor partial agonists. *J. Med. Chem.* **56**(10), 3852-3865 (2013).
2. Jones, R.W., Easter, R.A., McKeith, F.K., *et al.* Effect of the β-adrenergic agonist cimaterol (CL 263,780) on the growth and carcass characteristics of finishing swine. *J. Anim. Sci.* **61**(4), 905-913 (1985).
3. Byrem, T.M., Beermann, D.H., and Robinson, T.F. The β-agonist cimaterol directly enhances chronic protein accretion in skeletal muscle. *J. Anim. Sci.* **76**(4), 988-998 (1998).

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, 12/02/2022

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