

PRODUCT INFORMATION



Clencyclohexerol

Item No. 15365

CAS Registry No.: 157877-79-7

Formal Name: 4-amino-3,5-dichloro- α -[[[4-hydroxycyclohexyl)amino]methyl]-benzenemethanol

MF: $C_{14}H_{20}Cl_2N_2O_2$

FW: 319.2

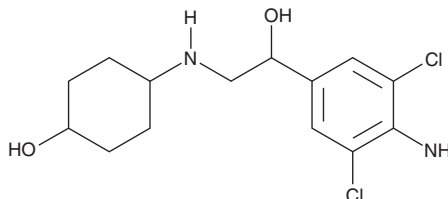
Purity: $\geq 90\%$

UV/Vis.: λ_{max} : 211, 248 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Clencyclohexerol is supplied as a crystalline solid. A stock solution may be made by dissolving the clencyclohexerol in the solvent of choice, which should be purged with an inert gas. Clencyclohexerol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of clencyclohexerol in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of clencyclohexerol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of clencyclohexerol in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Clencyclohexerol is an analog of clenbuterol (Item No. 14985). β -Adrenoceptor agonists, including clenbuterol, modulate protein synthesis and degradation and have therapeutic potential in muscle wasting disorders.¹ However, they are abused to stimulate muscle mass production, both in food-producing animals and in humans.^{2,3} Importantly, β -adrenoceptor agonists have deleterious cardiovascular side effects when used systemically and at high concentrations.¹ The physiological, cardiovascular, and toxicological properties of clencyclohexerol have not been evaluated. This product is intended for forensic and research purposes.

References

1. Ryall, J.G. and Lynch, G.S. The potential and the pitfalls of β -adrenoceptor agonists for the management of skeletal muscle wasting. *Pharmacol. Ther.* **120**(3), 219-232 (2008).
2. Salqu bre, G., Bresson, M., Villian, M., et al. Clenbuterol determination in calf hair by UPLC-MS-MS: Case report of a fraudulent use for cattle growth. *J. Anal. Toxicol.* **31**(2), 114-118 (2007).
3. Kanayama, G. and Pope, H.G., Jr. Illicit use of androgens and other hormones: Recent advances. *Curr. Opin. Endocrinol. Diabetes Obes.* **19**(3), 211-219 (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, 11/28/2022

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