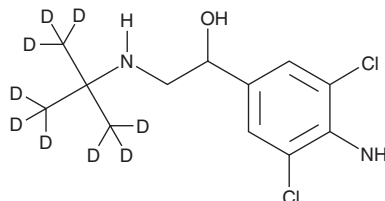


# PRODUCT INFORMATION



## Clenbuterol-d<sub>9</sub> Item No. 40816

**CAS Registry No.:** 129138-58-5  
**Formal Name:** 4-amino-3,5-dichloro-α-[[[1,1-di(methyl-d<sub>3</sub>)ethyl-2,2,2-d<sub>3</sub>]amino]methyl]-benzenemethanol  
**Synonym:** NAB 365-d<sub>9</sub>  
**MF:** C<sub>12</sub>H<sub>9</sub>Cl<sub>2</sub>D<sub>9</sub>N<sub>2</sub>O  
**FW:** 286.2  
**Chemical Purity:** ≥98% (clenbuterol)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>9</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Clenbuterol-d<sub>9</sub> is intended for use as an internal standard for the quantification of clenbuterol (Item No. 14985) 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).

Clenbuterol-d<sub>9</sub> is supplied as a solid. A stock solution may be made by dissolving the clenbuterol-d<sub>9</sub> in the solvent of choice, which should be purged with an inert gas. Clenbuterol-d<sub>9</sub> is slightly soluble in methanol and DMSO.

### Description

Clenbuterol is an agonist of β<sub>2</sub>-adrenergic receptors (β<sub>2</sub>-ARs).<sup>1</sup> It induces relaxation of equine tracheal muscle strips (EC<sub>50</sub> = 2.1 nM). Clenbuterol (300 nM) increases the number of autophagosomes and autophagic flux in HepG2 cells.<sup>2</sup> It increases blood flow to adipose tissue and induces weight gain in rats when administered at a dose of 2 mg/kg, as well as induces relaxation of bovine uterus.<sup>3,4</sup> Clenbuterol also inhibits infection of MDCK cells by an H1N1 influenza isolate *in vitro* (EC<sub>50</sub> = 9.4 μM).<sup>5</sup> Formulations containing clenbuterol have been used in the treatment of airway obstruction in horses.

### References

1. Törneke, K., Larsson, C.I., and Appelgren, L.E. A comparison between clenbuterol, salbutamol and terbutaline in relation to receptor binding and *in vitro* relaxation of equine tracheal muscle. *J. Vet. Pharmacol. Ther.* **21(5)**, 388-392 (1998).
2. Farah, B.L., Sinha, R.A., Wu, Y., *et al.* β-Adrenergic agonist and antagonist regulation of autophagy in HepG2 cells, primary mouse hepatocytes, and mouse liver. *PLoS One* **9(6)**, (2014).
3. Rothwell, N.J., Stock, M.J., and Sudera, D.K. Changes in tissue blood flow and β-receptor density of skeletal muscle in rats treated with the β<sub>2</sub>-adrenoceptor agonist clenbuterol. *Br. J. Pharmacol.* **90(3)**, 601-607 (1987).
4. Denooij, P.P. The use of clenbuterol for obstetrical procedures in forty cows and one horse. *Can. Vet. J.* **25(9)**, 357-359 (1984).
5. Jang, Y., Shin, J.S., Lee, J.-Y., *et al.* In vitro and in vivo antiviral activity of nylidrin by targeting the hemagglutinin 2-mediated membrane fusion of influenza A virus. *Viruses* **12(5)**, 581 (2020).

#### 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, 03/20/2024

#### 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