

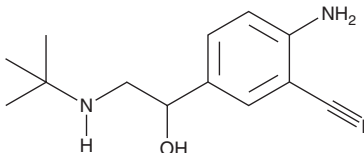
# PRODUCT INFORMATION



## Cimbuterol

Item No. 31743

**CAS Registry No.:** 54239-39-3  
**Formal Name:** 2-amino-5-[2-[(1,1-dimethylethyl)amino]-1-hydroxyethyl]-benzonitrile  
**MF:** C<sub>13</sub>H<sub>19</sub>N<sub>3</sub>O  
**FW:** 233.3  
**Purity:** ≥95%  
**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

Cimbuterol is supplied as a solid. A stock solution may be made by dissolving the cimbuterol in the solvent of choice, which should be purged with an inert gas. Cimbuterol is slightly soluble in methanol and dichloromethane.

### Description

Cimbuterol is an agonist of  $\beta$ -adrenergic receptors ( $\beta$ -ARs).<sup>1,2</sup> It has been used as a standard for the simultaneous detection of multiple veterinary residues in commercial animal meat by LC-MS/MS.<sup>1</sup> Cimbuterol has been used illicitly to promote growth and improve feed efficiency in livestock.

### References

1. Shao, B., Jia, X., Zhang, J., *et al.* Multi-residual analysis of 16  $\beta$ -agonists in pig liver, kidney and muscle by ultra performance liquid chromatography tandem mass spectrometry. *Food Chem.* **114(3)**, 1115-1121 (2009).
2. Robert, C., Gillard, N., Brasseur, P.-Y., *et al.* Rapid multi-residue and multi-class qualitative screening for veterinary drugs in foods of animal origin by UHPLC-MS/MS. *Food Addit. Contam. A.* **30(3)**, 443-457 (2013).

#### 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, 10/25/2022

#### 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