

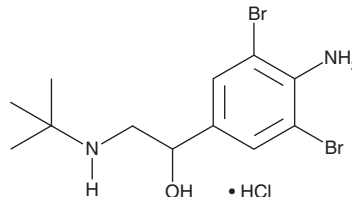
# PRODUCT INFORMATION



## Brombuterol (hydrochloride)

Item No. 33592

**CAS Registry No.:** 21912-49-2  
**Formal Name:** 4-amino-3,5-dibromo- $\alpha$ -[[[1,1-dimethylethyl]amino]methyl]-benzenemethanol, monohydrochloride  
**MF:** C<sub>12</sub>H<sub>18</sub>Br<sub>2</sub>N<sub>2</sub>O • HCl  
**FW:** 402.6  
**Purity:**  $\geq$ 95%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 215 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Brombuterol (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the brombuterol (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Brombuterol (hydrochloride) is soluble in DMSO and methanol.

### Description

Brombuterol is an agonist of the  $\beta_2$ -adrenergic receptor ( $\beta_2$ -AR).<sup>1</sup> It has been used as an illicit feed additive to increase the muscle-to-fat ratio of livestock.

### Reference

1. González-Antuña, A., Rodríguez-González, P., Centineo, G., *et al.* Simultaneous determination of seven  $\beta_2$ -agonists in human and bovine urine by isotope dilution liquid chromatography-tandem mass spectrometry using compound-specific minimally <sup>13</sup>C-labelled analogues. *J. Chromatogr. A.* **1372**, 63-71 (2014).

#### 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, 11/15/2022

#### 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