

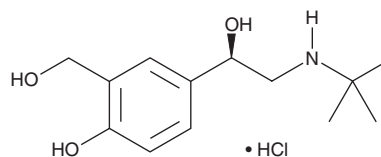
# PRODUCT INFORMATION



## (R)-Salbutamol (hydrochloride)

Item No. 23991

**CAS Registry No.:** 50293-90-8  
**Formal Name:**  $\alpha^1$ R-[[[1,1-dimethylethyl)amino]methyl]-4-hydroxy-1,3-benzenedimethanol, monohydrochloride  
**Synonyms:** Levalbuterol, (-)-Salbutamol,  
**MF:**  $C_{13}H_{21}NO_3 \cdot HCl$   
**FW:** 275.8  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 228, 276 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

(R)-Salbutamol (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the (R)-salbutamol (hydrochloride) in the solvent of choice. (R)-Salbutamol (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of (R)-salbutamol (hydrochloride) in these solvents is approximately 12, 20, and 25 mg/ml, respectively.

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 (R)-salbutamol (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of (R)-salbutamol (hydrochloride) in PBS, pH 7.2, is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

(R)-Salbutamol is an agonist of the  $\beta_2$ -adrenergic receptor ( $\beta_2$ -AR).<sup>1</sup> It induces relaxation of isolated guinea pig tracheal chains ( $EC_{50} = 16$  nM). (R)-Salbutamol selectively binds to the  $\beta_2$ -AR over the  $\beta_1$ -AR ( $K_{dS} = 0.236$  and  $1.54$   $\mu M$ , respectively).<sup>2</sup> Nebulized (R)-salbutamol reduces transpulmonary pressure in recurrent airway obstruction-affected horses ( $ED_{50} = 39.7$   $\mu g/animal$ ).<sup>3</sup> Formulations containing (R)-salbutamol have been used in the treatment or prevention of bronchospasm in patients with reversible obstructive airway disease.

### References

1. Hawkins, C.J. and Hawkins, G.T. Relative potency of (-)- and ( $\pm$ )-salbutamol on guinea pig tracheal tissue. *J. Med. Chem.* **16(7)**, 856-857 (1973).
2. Penn, R.B., Frielle, T., McCullough, J.R., *et al.* Comparison of R-, S-, and RS-albuterol interaction with human  $\beta_1$ - and  $\beta_2$ -adrenergic receptors. *Clin. Rev. Allergy Immunol.* **14(1)**, 37-45 (1996).
3. Arroyo, M.G., Cou  til, L.L., Nogradi, N., *et al.* Efficacy of inhaled levalbuterol compared to albuterol in horses with recurrent airway obstruction. *J. Vet. Intern. Med.* **30(4)**, 1333-1337 (2016).

#### 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, 03/03/2025

#### 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