

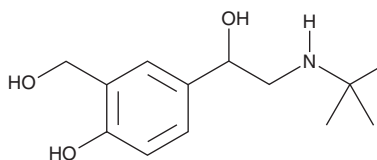
PRODUCT INFORMATION



Salbutamol

Item No. 21003

CAS Registry No.: 18559-94-9
Formal Name: α^1 -[[[1,1-dimethylethyl)amino)methyl]-4-hydroxy-1,3-benzenedimethanol
Synonyms: AH 3365, (\pm)-Salbutamol, DL-Salbutamol
MF: $C_{13}H_{21}NO_3$
FW: 239.3
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 226, 277 nm
Supplied as: A crystalline 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

Salbutamol is supplied as a crystalline solid. A stock solution may be made by dissolving the salbutamol in the solvent of choice. Salbutamol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of salbutamol in ethanol and DMF is approximately 10 mg/ml and approximately 5 mg/ml in DMSO.

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 salbutamol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of salbutamol in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Salbutamol is an agonist of the β_2 -adrenergic receptor (β_2 -AR).¹ It induces bronchodilation in isolated guinea pig trachea rings ($EC_{50} = 19$ nM). Salbutamol selectively binds to the β_2 -AR over the β_1 -AR and β_3 -AR in CHO cells expressing the human receptors (K_d s = 0.76, 21, and 46 μM , respectively).² Salbutamol (25 and 50 $\mu\text{g/kg}$, i.v.) reduces acetylcholine-induced bronchospasm in anesthetized guinea pigs.³ It also reduces the response of bronchial muscle to efferent vagal stimulation in anesthetized cats and dogs. Nebulized salbutamol reduces transpulmonary pressure in recurrent airway obstruction-affected horses ($sED_{50} = 43.6$ $\mu\text{g/animal}$).⁴ Formulations containing salbutamol have been used in the prevention of exercise-induced asthma and the prevention or treatment of chronic obstructive pulmonary disease (COPD).

References

1. Kern, C., Meyer, T., Droux, S., *et al.* Synthesis and pharmacological characterization of β_2 -adrenergic agonist enantiomers: Zilpaterol. *J. Med. Chem.* **52**(6), 1773-1777 (2009).
2. Baker, J.G. The selectivity of β -adrenoceptor antagonists at the human β_1 , β_2 and β_3 adrenoceptors. *Br. J. Pharmacol.* **144**(3), 317-322 (2005).
3. Cullum, V.A., Farmer, J.B., Jack, D., *et al.* Salbutamol: A new, selective β -adrenoceptive receptor stimulant. *Br. J. Pharmacol.* **35**(1), 141-151 (1969).
4. Arroyo, M.G., Couëtill, 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, 02/27/2025

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