

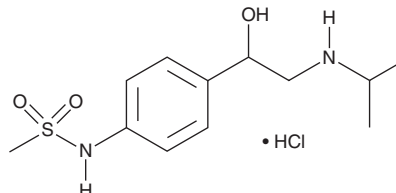
# PRODUCT INFORMATION



## Sotalol (hydrochloride)

Item No. 16136

**CAS Registry No.:** 959-24-0  
**Formal Name:** N-[4-[1-hydroxy-2-[(1-methylethyl)amino]ethyl]phenyl]-methanesulfonamide, monohydrochloride  
**Synonyms:** MJ1999, (±)-Sotalol, DL-Sotalol  
**MF:** C<sub>12</sub>H<sub>20</sub>N<sub>2</sub>O<sub>3</sub>S • HCl  
**FW:** 308.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 230 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

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

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 sotalol (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of sotalol (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Sotalol is a non-selective antagonist of  $\beta$ -adrenergic receptors ( $\beta$ -ARs; IC<sub>50</sub>s = 8.9 and 5.2  $\mu$ M for  $\beta$ 1- and  $\beta$ 2-ARs, respectively) and a class III antiarrhythmic agent.<sup>1,2</sup> It decreases delayed outward potassium currents (IK) in guinea pig ventricular cells and prolongs action potential duration in electrically stimulated isolated guinea pig papillary muscles when used at a concentration of 100  $\mu$ M.<sup>3</sup> Sotalol decreases heart rate and increases blood pressure and the cardiac functional refractory period (FRP) in a canine model of ventricular tachycardia induced by programmed electrical stimulation (PES).<sup>1</sup> Formulations containing sotalol have been used in the treatment of ventricular arrhythmias and maintenance of normal sinus rhythm in patients with atrial fibrillation or flutter (AFIB/AFL).

### References

1. Lis, R., Morgan, T.K., Jr., Marisca, A.J., et al. *J. Med. Chem.* **33**(10), 2883-2891 (1990).
2. Antonaccio, M.J. and Gomoll, A. *Am. J. Cardiol.* **65**(2), 12A-21A (1990).
3. Komeichi, K., Tohse, N., Nakaya, H., et al. *Eur. J. Pharm.* **187**(3), 313-322 (1990).

#### 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, 12/16/2022

#### 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