

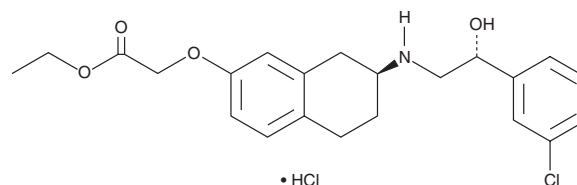
# PRODUCT INFORMATION



## SR 58611A (hydrochloride)

Item No. 11954

**CAS Registry No.:** 121524-09-2  
**Formal Name:** [[[7S)-7-[[[(2R)-2-(3-chlorophenyl)-2-hydroxyethyl]amino]-5,6,7,8-tetrahydro-2-naphthalenyl]oxy]acetic acid, ethyl ester, monohydrochloride  
**Synonym:** Amibegron  
**MF:** C<sub>22</sub>H<sub>26</sub>ClNO<sub>4</sub> • HCl  
**FW:** 440.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 277 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

SR 58611A (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 58611A (hydrochloride) in the solvent of choice. SR 58611A (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of SR 58611A (hydrochloride) in these solvents is approximately 30 mg/ml.

### Description

SR 58611A is a selective  $\beta_3$ -adrenergic receptor agonist ( $\beta_3$ -AR; EC<sub>50</sub> = 3.5 nM in rat colon).<sup>1</sup> Its activity cannot be blocked by the  $\beta_1$ - and  $\beta_2$ -AR antagonists CGP 20712A and ICI 118551 (Item No. 15591), respectively. SR 58611A is minimally active against  $\beta_1$ -ARs in rat uterus (EC<sub>50</sub> = 499 nM) and inactive against  $\beta_2$ -ARs in guinea pig trachea and atrium (EC<sub>50</sub>s = >10,000 and >30,000 nM, respectively). SR 58611A activates brown fat metabolism through activation of adenylyl cyclase activity and glycerol release in brown adipocytes (EC<sub>50</sub>s = 20 and 11 nM, respectively).<sup>2</sup> It also reduces hypothermia produced by apomorphine (Item No. 16094) and reserpine (Item No. 16474) and potentiates toxicity produced by yohimbine (Item No. 19869) in mice.<sup>3</sup> SR 58611A (0.6 to 2 mg/kg/day) also reduces the number of escape failures in a learned helplessness model of antidepressant-like activity in rats without changes in locomotor activity typically seen with  $\beta_2$ -AR agonists.

### References

1. Bianchetti, A. and Manara, L. *In vitro* inhibition of intestinal motility by phenylethanolaminotetralines: Evidence of atypical  $\beta$ -adrenoceptors in rat colon. *Br. J. Pharmacol.* **100**(4), 831-839 (1990).
2. Nisoli, E., Tonello, C., and Carruba, M.O. SR 58611A: A novel thermogenic  $\beta$ -adrenoceptor agonist. *Eur. J. Pharmacol.* **259**(2), 181-186 (1994).
3. Simiand, J., Keane, P.E., Guitard, J., *et al.* Antidepressant profile in rodents of SR 58611A, a new selective agonist for atypical  $\beta$ -adrenoceptors. *Eur. J. Pharmacol.* **219**(2), 193-201 (1992).

#### 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, 08/08/2017

#### 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