

# PRODUCT INFORMATION

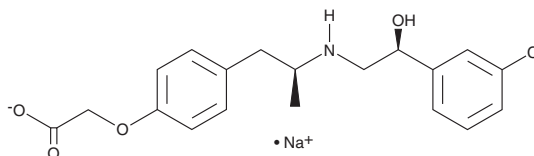


## BRL 37344 (sodium salt)

Item No. 18622

**CAS Registry No.:** 127299-93-8  
**Formal Name:** *rel*-2-[4-[(2*S*)-2-[[[(2*S*)-2-(3-chlorophenyl)-2-hydroxyethyl]amino]propyl]phenoxy]-acetic acid, monosodium salt

**Synonym:** BRL 37344A  
**MF:** C<sub>19</sub>H<sub>21</sub>ClNO<sub>4</sub> • Na  
**FW:** 385.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 274 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

BRL 37344 (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the BRL 37344 (sodium salt) in the solvent of choice, which should be purged with an inert gas. BRL 37344 (sodium salt) is soluble in the organic solvent DMSO at a concentration of approximately 5 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 BRL 37344 (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of BRL 37344 (sodium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

BRL 37344 is a selective agonist of  $\beta_3$ -adrenergic receptors with K<sub>i</sub> values of 0.43, 9.17, and 37.9  $\mu$ M for  $\beta_3$ ,  $\beta_2$ , and  $\beta_1$ , respectively.<sup>1</sup>  $\beta_3$ -Adrenoceptors are expressed in adipose tissue as well as endothelial and myocardial tissues.<sup>2</sup> This compound has been used to explore the role of these receptors in the regulation of thermogenesis and lipolysis and to modulate cardiac function in models of cardiovascular disease.<sup>3</sup> It has also been used in rodents to study the role that  $\beta_3$ -adrenergic receptors play in regulating anxiety-like behaviors.<sup>4</sup>

### References

- Hoffmann, C., Leitz, M.R., Oberdorf-Maass, S., et al. Comparative pharmacology of human  $\beta$ -adrenergic receptor subtypes-characterization of stably transfected receptors in CHO cells. *Naunyn Schmiedeberg's Arch. Pharmacol.* **369**(2), 151-159 (2004).
- Kanzler, S.A., Januario, A.C., and Paschoalini, M.A. Involvement of  $\beta_3$ -adrenergic receptors in the control of food intake in rats. *Braz. J. Med. Biol. Res.* **44**(11), 1141-1147 (2011).
- Niu, X., Zhao, L., Li, X., et al.  $\beta_3$ -Adrenoreceptor stimulation protects against myocardial infarction injury via eNOS and nNOS activation. *PLoS One* **9**(6), (2014).
- Butler, T.R., Chappell, A.M., and Weiner, J.L. Effect of  $\beta_3$  adrenoceptor activation in the basolateral amygdala on ethanol seeking behaviors. *Psychopharmacology (Berl)* **231**(1), 293-303 (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, 12/14/2022

#### 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