

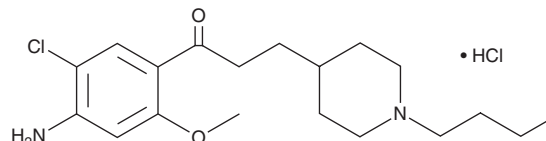
PRODUCT INFORMATION



RS 67333 (hydrochloride)

Item No. 45937

CAS Registry No.: 168986-60-5
Formal Name: 1-(4-amino-5-chloro-2-methoxyphenyl)-3-(1-butyl-4-piperidinyl)-1-propanone, monohydrochloride
MF: $C_{19}H_{29}ClN_2O_2 \cdot HCl$
FW: 389.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

RS 67333 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the RS 67333 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RS 67333 (hydrochloride) is soluble (≥ 10 mg/ml) in DMSO.

Description

RS 67333 is an agonist of the serotonin (5-HT) receptor subtype 5-HT₄.¹ It selectively binds to 5-HT₄ ($K_i = 2$ nM) over 5-HT_{1A}, 5-HT_{2A}, 5-HT_{2C}, and 5-HT₃ receptors (K_i s = 398, 501, 794, and 398 nM, respectively), as well as α - and β -adrenergic and muscarinic receptors (K_i s = >200 nM for all). It also binds to σ_1 - and σ_2 -receptors (K_i s = 1 and 10 nM, respectively). RS 67333 induces relaxation of carbachol-precontracted rat esophageal muscularis mucosae preparations ($EC_{50} = 2$ nM) and induces tachycardia in anesthetized micropigs at 4.9 μ g/kg. It reverses scopolamine-induced memory deficits in mice in the Y-maze test when administered at a dose of 1 mg/kg.²

References

1. Eglén, R.M., Bonhaus, D.W., Johnson, L.G., *et al.* Pharmacological characterization of two novel and potent 5-HT₄ receptor agonists, RS 67333 and RS 67506, *in vitro* and *in vivo*. *Br. J. Pharmacol.* **115**(8), 1387-1392 (1995).
2. Lelong, V., Lhonnear, L., Dauphin, F., *et al.* BIMU 1 and RS 67333, two 5-HT₄ receptor agonists, modulate spontaneous alternation deficits induced by scopolamine in the mouse. *Naunyn Schmiedeberg's Arch Pharmacol.* **367**(6), 621-628 (2003).

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, 07/02/2026

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