

PRODUCT INFORMATION

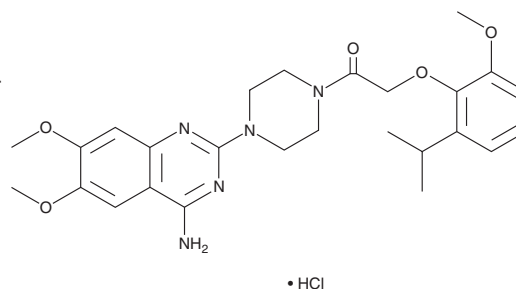


Rec 15/2615 (hydrochloride)

Item No. 27677

CAS Registry No.: 173059-17-1
Formal Name: 1-[4-(4-amino-6,7-dimethoxy-2-quinazoliny)-1-piperaziny]-2-[2-methoxy-6-(1-methylethyl)phenoxy]-ethanone, monohydrochloride

Synonym: SB 216469
MF: C₂₆H₃₃N₅O₅ • HCl
FW: 532.0
Purity: ≥98%
Supplied as: A 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

Rec 15/2615 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the rec 15/2615 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Rec 15/2615 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

Rec 15/2615 is an antagonist of α_{1B} -adrenergic receptors (α_{1B} -ARs; $K_i = 0.45$ nM for the recombinant human receptor).¹ It selectively inhibits α_{1B} -ARs over α_{1A} -, α_{1D} -, and α_{1L} -ARs (K_i s = 7.59, 10.23, and 49 nM, respectively). Rec 15/2615 inhibits norepinephrine-induced contractions of isolated rabbit prostate and urethral strips (K_i s = 100 and 316.2 nM, respectively), as well as reduces norepinephrine-induced contractions of chloroethylclonidine-precontracted isolated rabbit aortic rings ($K_i = 50$ nM).² It decreases diastolic blood pressure ($ED_{25} = 183$ μ g/kg, i.v.) and increases intracavernous pressure in anesthetized dogs when administered intracavernously at doses ranging from 30 and 1,000 μ g/kg.^{1,2}

References

1. Sironi, G., Colombo, D., Poggesi, E., *et al.* Effects of intracavernous administration of selective antagonists of α_1 -adrenoceptor subtypes on erection in anesthetized rats and dogs. *J. Pharmacol. Exp. Ther.* **292**(3), 974-981 (2000).
2. Testa, R., Guarneri, L., Angelico, P., *et al.* Pharmacological characterization of the uroselective *alpha*-1 antagonist Rec 15/2739 (SB 216469): Role of the *alpha*-1L adrenoceptor in tissue selectivity, part II. *J. Pharmacol. Exp. Ther.* **281**(3), 1284-1293 (1997).

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, 03/05/2026

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