

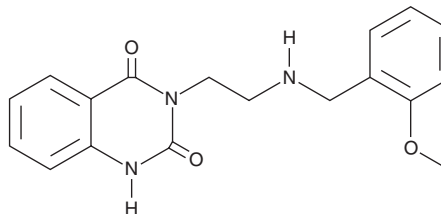
PRODUCT INFORMATION



RH-34

Item No. 14445

CAS Registry No.: 1028307-48-3
Formal Name: 3-[2-[[[(2-methoxyphenyl)methyl]amino]ethyl]-2,4(1H,3H)-quinazolidione
MF: C₁₈H₁₉N₃O₃
FW: 325.4
Purity: ≥95%
UV/Vis.: λ_{max}: 219, 278, 312 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Ketanserin is a potent antagonist of the serotonin receptor 5-HT_{2A} (K_i = 3.16 nM) that has hypertensive effects *in vivo*.^{1,2} RH-34 is a derivative of ketanserin in which the 4-(*p*-fluorobenzoyl)piperidine group has been replaced with an N-(2-methoxybenzyl) moiety. In phenethylamines, which are 5-HT_{2A} agonists, this addition increases the affinity and selectivity for the 5-HT_{2A} receptor over other serotonin receptors.³ The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

References

1. Bonhaus, D.W., Bach, C., DeSouza, A., *et al.* The pharmacology and distribution of human 5-hydroxytryptamine_{2B} (5-HT_{2B}) receptor gene products: Comparison with 5-HT_{2A} and 5-HT_{2C} receptors. *Br. J. Pharmacol.* **115(4)**, 622-628 (1995).
2. van Zwieten, P.A., Blauw, G.J., and van Brummelen, P. The role of 5-hydroxytryptamine and 5-hydroxytryptaminergic mechanisms in hypertension. *Br. J. Clin. Pharmacol.* **30(Suppl 1)**, 69S-74S (1990).
3. Braden, M.R., Parrish, J.C., Naylor, J.C., *et al.* Molecular interaction of serotonin 5-HT_{2A} receptor residues Phe339^{6.51} and Phe340^{6.52} with superpotent N-benzyl phenethylamine agonists. *Mol. Pharm.* **70(6)**, 1956-1965 (2006).

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, 06/19/2023

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