

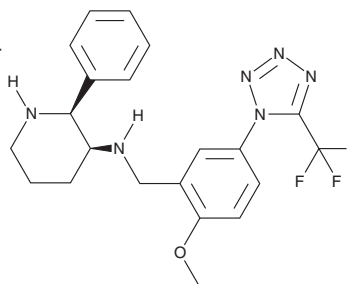
PRODUCT INFORMATION



Vofopitant

Item No. 42029

CAS Registry No.: 168266-90-8
Formal Name: N-[[2S-methoxy-5-[5-(trifluoromethyl)-1H-tetrazol-1-yl]phenyl]methyl]-2-phenyl-3S-piperidinamine
Synonym: GR 205171
MF: C₂₁H₂₃F₃N₆O
FW: 432.4
Purity: ≥95%
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

Vofopitant is supplied as a solid. A stock solution may be made by dissolving the vofopitant in the solvent of choice, which should be purged with an inert gas. Vofopitant is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Vofopitant is a neurokinin-1 (NK₁) receptor antagonist (K_i = 0.03 nM for the human receptor).¹ It is selective for NK₁ over NK₂ and NK₃ receptors (IC₅₀s = >10,000 nM for both), the serotonin (5-HT) receptor subtypes 5-HT_{1A}, 5-HT_{1D}, and 5-HT_{2A} (K_is = 501, 251, and 316 nM, respectively), and histamine H₁ and H₂ receptors (K_is = 316 and 251 nM, respectively). It inhibits systemic vascular leakage induced by resiniferatoxin in guinea pigs (ID₅₀ = 0.007 mg/kg).² Vofopitant (0.1 mg/kg) inhibits retching and vomiting induced by various emetogens in ferrets.¹ It increases the number of entries into the open arms of the elevated plus maze in gerbils when administered at doses ranging from 0.3 to 5 mg/kg and increases the percentage of time spent in the open arms at 5 mg/kg, indicating anxiolytic-like activity.³ Vofopitant potentiates the reduction in immobility time induced by citalopram (Item No. 14572) in the forced swim test in mice when administered at a dose of 40 mg/kg but inhibits the reduction in social interaction induced by citalopram in rats and gerbils at 0.63 and 10 mg/kg, indicating potentiation and inhibition of citalopram-induced antidepressant-like and anxiety-like behaviors, respectively.⁴

References

1. Gardner, C.J., Armour, D.R., Beattie, D.T., *et al.* GR205171: A novel antagonist with high affinity for the tachykinin NK₁ receptor, and potent broad-spectrum anti-emetic activity. *Regul. Pept.* **65(1)**, 45-53 (1996).
2. Hale, J.J., Mills, S.G., MacCoss, M., *et al.* Structural optimization affording 2-(R)-(1-(R)-3, 5-bis (trifluoromethyl)phenylethoxy)-3-(S)-(4-fluoro)phenyl-4-(3-oxo-1,2,4-triazol-5-yl)methylmorpholine, a potent, orally active, long-acting morpholine acetal human NK-1 receptor antagonist. *J. Med. Chem.* **41(23)**, 4607-4614 (1998).
3. Heldt, S.A., Davis, M., Ratti, E., *et al.* Anxiolytic-like effects of the neurokinin 1 receptor antagonist GR-205171 in the elevated plus maze and contextual fear-potentiated startle model of anxiety in gerbils. *Behav. Pharmacol.* **20(7)**, 584-595 (2009).
4. Gobert, A., Brocco, M., Dekeyne, A., *et al.* Neurokinin₁ antagonists potentiate antidepressant properties of serotonin reuptake inhibitors, yet blunt their anxiogenic actions: A neurochemical, electrophysiological, and behavioral characterization. *Neuropsychopharmacology* **34(4)**, 1039-1056 (2009).

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, 10/25/2024

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