

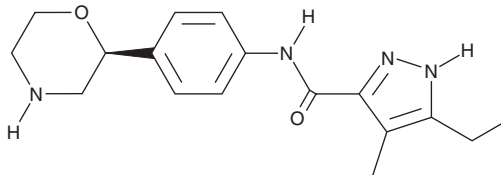
PRODUCT INFORMATION



Ralmitaront

Item No. 41111

CAS Registry No.: 2133417-13-5
Formal Name: 5-ethyl-4-methyl-N-[4-(2S)-2-morpholinylphenyl]-1H-pyrazole-3-carboxamide
Synonyms: RG-7906, RO6889450
MF: C₁₇H₂₂N₄O₂
FW: 314.4
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

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

Description

Ralmitaront is an agonist of trace amine-associated receptor 1 (TAAR1).¹ It induces cAMP accumulation in HEK293T cells expressing human TAAR1 (EC₅₀ = 17.78 nM).

Reference

1. Ågren, R., Betari, N., Saarinen, M., *et al.* In vitro comparison of ulotaront (SEP-363856) and ralmitaront (RO6889450): Two TAAR1 agonist candidate antipsychotics. *Int. J. Neuropsychopharmacol.* **26(9)**, 599-606 (2023).

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, 06/13/2024

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