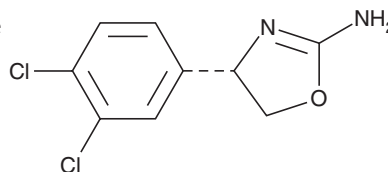


PRODUCT INFORMATION



RO5203648
Item No. 45362

CAS Registry No.: 1043491-54-8
Formal Name: 4-(3,4-dichlorophenyl)-4,5-dihydro-2-oxazolamine
MF: C₉H₈Cl₂N₂O
FW: 231.1
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

RO5203648 is supplied as a solid. A stock solution may be made by dissolving the RO5203648 in the solvent of choice, which should be purged with an inert gas. RO5203648 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

RO5203648 is a partial agonist of trace amine-associated receptor 1 (TAAR1).¹ It increases cAMP production in HEK293 cells expressing TAAR1 (EC₅₀s = 4, 6.8, 30, and 31 nM for the mouse, rat, human, and monkey receptors, respectively). RO5203648 (500 nM) increases the firing frequency of dopaminergic and serotonergic neurons in the ventral tegmental area (VTA) and dorsal raphe nucleus (DRN) in brain slices from wild-type but not *Taar1*^{-/-} mice. It reduces cocaine-induced hyperlocomotion in mice when administered at doses of 1 and 3, but not 10, mg/kg. It also reduces catalepsy induced by haloperidol (Item No. 12014) and reduces immobility time in the forced swim test in rats. RO5203648 (10 mg/kg) increases the percentage of correct responses in an object retrieval task in cynomolgus monkeys and reduces cocaine administration in a self-administration paradigm in rats.

Reference

1. Revel, F.G., Moreau, J.L., Gainetdinov, R.R., *et al.* Trace amine-associated receptor 1 partial agonism reveals novel paradigm for neuropsychiatric therapeutics. *Biol. Psychiatry* **72(11)**, 934-942 (2012).

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, 04/21/2026

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