

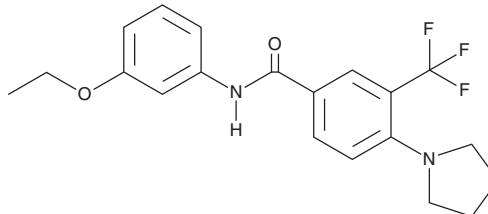
PRODUCT INFORMATION



EPPTB

Item No. 34330

CAS Registry No.: 1110781-88-8
Formal Name: N-(3-ethoxyphenyl)-4-(1-pyrrolidinyl)-3-(trifluoromethyl)-benzamide
Synonym: Ro 5212773
MF: C₂₀H₂₁F₃N₂O₂
FW: 378.4
Purity: ≥90%
UV/Vis.: λ_{max}: 290 nm
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

EPPTB is supplied as a solid. A stock solution may be made by dissolving the EPPTB in the solvent of choice, which should be purged with an inert gas. EPPTB is soluble in ethanol.

Description

EPPTB is an antagonist of trace amine-associated receptor 1 (TAAR1; K_i = 0.9 nM for the mouse receptor).¹ It selectively inhibits β-phenylethylamine-induced cAMP production in HEK293 cells expressing mouse TAAR1 over human or rat TAAR1 (IC₅₀s = 27.5, 7,487, and 4,539 nM, respectively). EPPTB (10 nM) reduces decreases in the firing frequency of dopaminergic neurons induced by *p*-tyramine (Item No. 18601) in mouse ventral tegmental area (VTA) slices when used at a concentration of 10 nM.

Reference

1. Bradaia, A., Trube, G., Stalder, H., *et al.* The selective antagonist EPPTB reveals TAAR1-mediated regulatory mechanisms in dopaminergic neurons of the mesolimbic system. *Proc. Natl. Acad. Sci. USA* **106**(47), 20081-20086 (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, 05/09/2023

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