

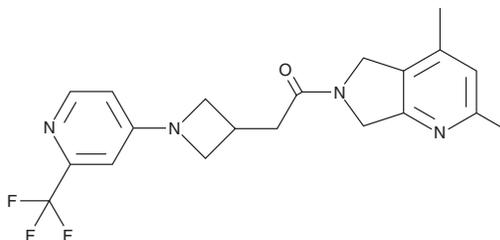
PRODUCT INFORMATION



Emraclidine

Item No. 42125

CAS Registry No.: 2170722-84-4
Formal Name: 1-(5,7-dihydro-2,4-dimethyl-6H-pyrrolo[3,4-b]pyridin-6-yl)-2-[1-[2-(trifluoromethyl)-4-pyridinyl]-3-azetidiny]-ethanone
Synonyms: CVL-231, PF-06852231
MF: C₂₀H₂₁F₃N₄O
FW: 390.4
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

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

Description

Emraclidine is a positive allosteric modulator (PAM) of M₄ muscarinic acetylcholine receptors (mAChRs).¹ It is selective for M₄ mAChRs over a panel of 59 receptors, ion channels, transporters, and enzymes at 10 μM. Emraclidine selectively potentiates ACh-induced activation of the M₄ mAChR over M₁-M₃ and M₅ mAChRs in cells expressing the human receptors (EC₅₀s = 12.3, 4,798, >10,000, >10,000, and >10,000 nM, respectively). *In vivo*, emraclidine (3.2 mg/kg) inhibits amphetamine-induced locomotor activity in mice and prepulse inhibition deficits in rats.

Reference

1. Butler, C.R., Popiolek, M., McAllister, L.A., *et al.* Design and synthesis of clinical candidate PF-06852231 (CVL-231): A brain penetrant, selective, positive allosteric modulator of the M4 muscarinic acetylcholine receptor. *J. Med. Chem.* **67**(13), 10831-10847 (2024).

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, 10/29/2024

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