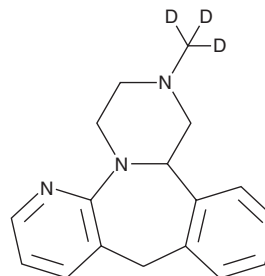


PRODUCT INFORMATION



Mirtazapine-d₃ Item No. 21126

CAS Registry No.: 1216678-68-0
Formal Name: 1,2,3,4,10,14b-hexahydro-2-methyl-d₃-pyrazino[2,1-a]pyrido[2,3-c][2]benzazepine
Synonym: 6-Azamianserin-d₃
MF: C₁₇H₁₆D₃N₃
FW: 268.4
Chemical Purity: ≥98% (Mirtazapine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A neat solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Mirtazapine-d₃ (Item No. 21126) is an analytical reference material intended for use as an internal standard for the quantification of mirtazapine (Item No. 19994). The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Mirtazapine is categorized as an antidepressant.¹ It inhibits maintenance of methamphetamine-induced conditioned place preference in rats. It also impairs acquisition and reinstatement of cocaine-induced place preference in rats.² This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Voigt, R.M., Mickiewicz, A.L., Napier, T.C. Repeated mirtazapine nullifies the maintenance of previously established methamphetamine-induced conditioned place preference in rats. *Behav. Brain Res.* **225(1)**, 91-96 (2011).
2. Barbosa-Méndez, S., Matus-Ortega, M., Jacinto-Gutiérrez, S., et al. Mirtazapine impairs acquisition and reinstatement of cocaine-induced place preference in rats. *Eur. J. Pharmacol.* **820**, 183-190 (2018).

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, 08/21/2020

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