

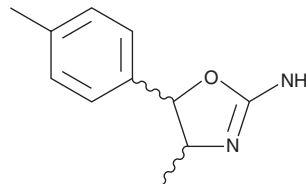
PRODUCT INFORMATION



4,4'-Dimethylaminorex

Item No. 15281

CAS Registry No.: 1445569-01-6
Formal Name: 4,5-dihydro-4-methyl-5-(4-methylphenyl)-2-oxazolamine
Synonyms: 4,4'-DMAR, 4,4'-DMAX
MF: C₁₁H₁₄N₂O
FW: 190.3
Purity: ≥98% (mixture of diastereomers)
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

4,4'-Dimethylaminorex (4,4'-DMAR) is a designer drug and *para*-methyl derivative of 4-methylaminorex, an amphetamine-like stimulant.^{1,2} 4,4'-DMAR is a less common alternative to 4-methylaminorex, although the latter is an internationally controlled substance.³ The physiological and toxicological properties of this compound are not known. 4,4'-DMAR is regulated as a Schedule I compound in the United States. This product is intended for forensic and research applications.

References

1. Hanson, G.R., Bunker, C.F., Johnson, M., *et al.* Response of monoaminergic and neuropeptide systems to 4-methylaminorex: A new stimulant of abuse. *Eur. J. Pharmacol.* **218(2-3)**, 287-293 (1992).
2. Kankaanpää, A., Ellermaa, S., Meririnne, E., *et al.* Acute neurochemical and behavioral effects of stereoisomers of 4-methylaminorex in relation to brain drug concentrations. *J. Pharmacol. Exp. Ther.* **300(2)**, 450-459 (2002).
3. Nizar, H., Dargan, P.I., and Wood, D.M. Using internet snapshot surveys to enhance our understanding of the availability of the novel psychoactive substance 4-methylaminorex and 4,4'-dimethylaminorex. *J. Med. Toxicol.* **11(1)**, 80-84 (2014).

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, 03/29/2024

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