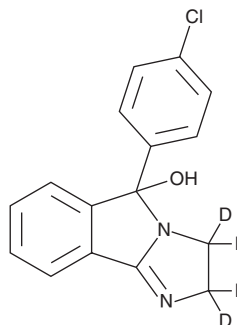


PRODUCT INFORMATION



Mazindol-d₄ Item No. 21640

CAS Registry No.: 1246815-50-8
Formal Name: 5-(4-chlorophenyl)-2,5-dihydro-3H-imidazo[2,1-a]isoindol-5-ol-d₄
MF: C₁₆H₉ClD₄N₂O
FW: 288.8
Chemical Purity: ≥98% (Mazindol)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
UV/Vis.: λ_{max}: 225, 268 nm
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

Mazindol-d₄ (Item No. 21640) is intended for use as an internal standard for the quantification of mazindol (Item No. 21642) by GC- or LC-MS. 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).

Mazindol is categorized as a stimulant and can be detected in urine and plasma.¹ It blocks reuptake at the norepinephrine and dopamine transporters, with less inhibition at the serotonin transporter (K_s = 3.2, 27.6, and 153 nM, respectively).² Mazindol is used illegally as an anorectic. This product is intended for research and forensic applications.

References

1. de Oliveira, M.H., Ferreira, P.C.L., Carlos, G., *et al.* Validation and application of a liquid chromatography-electrospray ionization mass spectrometric method for determination of mazindol in human plasma and urine. *J. Pharmacol. Toxicol. Methods* **79**, 1-6 (2016).
2. Zhou, J. Norepinephrine transporter inhibitors and their therapeutic potential. *Drugs Future* **29(12)**, 1235-1244 (2004).

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, 07/06/2023

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