

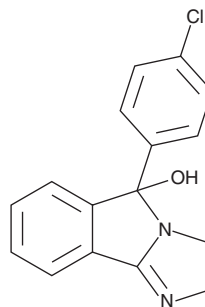
PRODUCT INFORMATION



Mazindol

Item No. 21642

CAS Registry No.: 22232-71-9
Formal Name: 5-(4-chlorophenyl)-2,5-dihydro-3H-imidazo[2,1-a]isoindol-5-ol
Synonyms: AN 448, SAH 42548
MF: C₁₆H₁₃ClN₂O
FW: 284.7
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 266 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 (Item No. 21642) is an analytical reference standard structurally categorized as a stimulant that 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 Formulations containing mazindol were previously used in treatment of Duchenne muscular dystrophy but it was discontinued from use in the United States. 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