

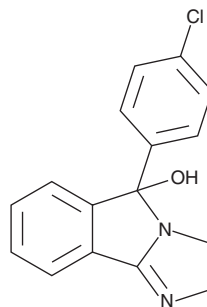
PRODUCT INFORMATION



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 categorized as a stimulant and inhibitor of norepinephrine, dopamine, and serotonin reuptake.¹ Mazindol decreases high-fat diet-induced weight gain in mice without eliciting abuse potential or increasing locomotor activity.² Formulations containing mazindol have previously been used in the management of obesity. Mazindol is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

References

1. Konofal, E., Arnulf, I., Bizot, J.C., *et al.* Mazindol immediate-release/sustained-release (IR/SR): A 50-Year legacy of multifaceted mechanisms and emerging therapeutic potential. *Clin. Drug Investig.* **46**(3), 259-280 (2026).
2. Aotani, D., Son, C., Shimizu, Y., *et al.* Reevaluation of anti-obesity action of mazindol and elucidation of its effect on the reward system. *Neurosci. Lett.* **633**, 141-145 (2016).

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, 06/30/2026

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