

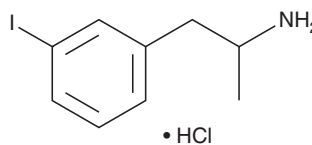
PRODUCT INFORMATION



3-Iodoamphetamine (hydrochloride)

Item No. 9001852

CAS Registry No.: 20110-56-9
Formal Name: 3-iodo- α -methyl-benzeneethanamine, monohydrochloride
Synonyms: *m*-iodo- α -methyl-Phenethylamine, 3-IA, *m*-IA
MF: C₉H₁₂I_N • HCl
FW: 297.6
Purity: ≥98%
UV/Vis.: λ_{max} : 231 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

Some halogenated amphetamines deplete brain serotonin levels and display pronounced neurotoxicity.¹ 3-Iodoamphetamine is a halogenated amphetamine featuring an iodine atom at the *meta* position of the phenyl group. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

Reference

1. Fuller, R.W., Baker, J.C., Perry, K.W., *et al.* Comparison of 4-chloro-, 4-bromo- and 4-fluoroamphetamine in rats: Drug levels in brain and effects on brain serotonin metabolism. *Neuropharmacology* **14**(10), 739-746 (1975).

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, 07/12/2023

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