

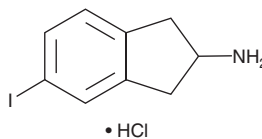
PRODUCT INFORMATION



5-IAI (hydrochloride)

Item No. 11035

CAS Registry No.: 1782044-60-3
Formal Name: 2,3-dihydro-5-iodo-1H-inden-2-amine, monohydrochloride
Synonym: 5-iodo-2-Aminoindan
MF: C₉H₁₀IN • HCl
FW: 295.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Aminoindans (AI) were originally studied as semi-rigid congeners of phenylethylamines, which are psychoactive alkaloids.^{1,2} 5-IAI is a psychoactive analog of *p*-iodoamphetamine and is indistinguishable, in its physiological effects, from 3,4-methylenedioxymethamphetamine in rats.³ A single 40 mg/kg dose of 5-IAI significantly reduces both serotonin uptake sites and hippocampal serotonin levels in rats.⁴ This product is intended for forensic applications.

References

1. Cannon, J.G., Perez, J.A., and Pease, J.P. Comparison of biological effects of N-alkylated congeners of β -phenethylamine derived from 2-aminotetralin, 2-aminoindan, and 6-aminobenzocycloheptene. *J. Med. Chem.* **23**, 745-749 (1980).
2. Nichols, D.E., Brewster, W.K., Johnson, M.P., et al. Nonneurotoxic tetralin and indan analogues of 3,4-(methylenedioxy)amphetamine (MDA). *J. Med. Chem.* **33(2)**, 703-710 (1990).
3. Johnson, M.P., Conarty, P.F., and Nichols, D.E. [3H]monoamine releasing and uptake inhibition properties of 3,4-methylenedioxymethamphetamine and *p*-chloroamphetamine analogues. *Eur. J. Pharmacol.* **200**, 9-16 (1991).
4. Nichols, D.E., Johnson, M.P., and Oberlender, R. 5-Iodo-2-aminoindan, a nonneurotoxic analogue of *p*-iodoamphetamine. *Pharmacol. Biochem. Behav.* **38**, 135-139 (1991).

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, 11/30/2017

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