

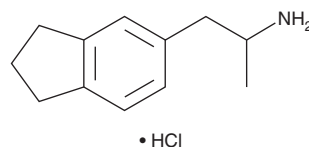
PRODUCT INFORMATION



5-APDI (hydrochloride)

Item No. 13899

CAS Registry No.: 152623-95-5
Formal Name: 2,3-dihydro- α -methyl-1H-indene-5-ethanamine, monohydrochloride
Synonyms: 5-(2-Aminopropyl)-2,3-dihydro-1H-indene, IAP, Indanylamino propane
MF: C₁₂H₁₇N • HCl
FW: 211.7
Purity: ≥98%
UV/Vis.: λ_{max} : 271, 277 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

5-APDI is an indane with an amphetamine-like aminopropane group. It inhibits the uptake of serotonin, dopamine, and norepinephrine by crude synaptosomes (IC_{50} = 82, 1,847, and 849 nM, respectively).¹ In discrimination studies using rats, 5-APDI fully substitutes for the entactogen N-methyl-1,3-dioxolyl-N-methylbutanamine but does not substitute for amphetamine.¹ This product is intended for forensic and research applications.

Reference

1. Monte, A.P., Marona-Lewicka, D., Cozzi, N.V., *et al.* Synthesis and pharmacological examination of benzofuran, indan, and tetralin analogues of 3,4-(methylenedioxy)amphetamine. *J. Med. Chem.* **36**(23), 3700-3706 (1993).

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, 07/25/2023

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