

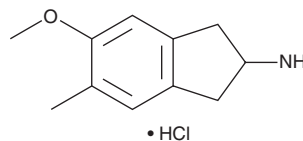
PRODUCT INFORMATION



MMAI (hydrochloride)

Item No. 13895

CAS Registry No.: 132980-17-7
Formal Name: 2,3-dihydro-5-methoxy-6-methyl-1H-inden-2-amine, monohydrochloride
MF: $C_{11}H_{15}NO \cdot HCl$
FW: 213.7
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 284 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

MMAI is a psychoactive aminoindan which can be found in recreational drug mixtures.¹ It potently inhibits serotonin uptake ($IC_{50} = 212$ nM) and stimulates its release without inducing the neurotoxicity induced by 3,4-methylenedioxy methamphetamine and similar compounds.^{2,3} This product is intended for forensic and research applications.

References

1. Sainsbury, P.D., Kicman, A.T., Archer, R.P., *et al.* Aminoindanes--the next wave of 'legal highs'? *Drug Test Anal.* **3**(7-8), 479-482 (2011).
2. Johnson, M.P., Frescas, S.P., Oberlender, R., *et al.* Synthesis and pharmacological examination of 1-(3-methoxy-4-methylphenyl)-2-aminopropane and 5-methoxy-6-methyl-2-aminoindan: Similarities to 3,4-(methylenedioxy)methamphetamine (MDMA). *J. Med. Chem.* **34**(5), 1662-1668 (1991).
3. 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