# **PRODUCT** INFORMATION



**MEAI** (hydrochloride)

Item No. 18286

CAS Registry No.:	81593-54-6
Formal Name:	2,3-dihydro-5-methoxy-1H-Inden-2-amine,
	monohydrochloride
Synonyms:	5-methoxy-2-Al,
	5-methoxy-2-Aminoindane,
	5-methoxy Al
MF:	C <sub>10</sub> H <sub>13</sub> NO • HCl
FW:	199.7
Purity:	≥95%
UV/Vis.:	λ <sub>max</sub> : 225, 281 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

Aminoindanes (Als) were originally studied as semi-rigid congeners of phenylethylamines, which are psychoactive alkaloids.<sup>1-3</sup> MEAI is a derivative of MMAI (Item No. 13895), a psychoactive aminoindane that has been identified in recreational drug mixtures.<sup>1</sup> MEAI differs in structure from MMAI by the absence of a methyl group. This product is intended for forensic and research purposes.

## 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. 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(7), 745-749 (1980).
- 3. 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).

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.

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, 09/20/2023

## 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