

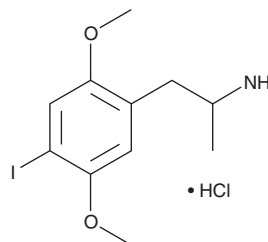
PRODUCT INFORMATION



DOI (hydrochloride)

Item No. 13885

CAS Registry No.: 42203-78-1
Formal Name: 4-iodo-2,5-dimethoxy- α -methylbenzeneethanamine, monohydrochloride
Synonym: 2,5-dimethoxy-4-Iodoamphetamine
MF: C₁₁H₁₆INO₂ • HCl
FW: 357.6
Purity: ≥98%
UV/Vis.: λ_{max} : 235, 300 nm
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

DOI (hydrochloride) (Item No. 13885) is an analytical reference standard categorized as an amphetamine and phenethylamine. DOI has psychotomimetic properties and abuse of DOI has been reported.¹⁻³ This product is intended for research and forensic applications.

References

1. Shulgin, A.T. Psychotomimetic Drugs: Structure-activity relationships. *Stimulants* 243-333 (1978).
2. Bruno, R., Matthews, A.J., Dunn, M., *et al.* Emerging psychoactive substance use among regular ecstasy users in Australia. *Drug Alcohol Depend.* **124**(1-2), 19-25 (2012).
3. Kerrigan, S., Banuelos, S., Perrella, L., *et al.* Simultaneous detection of ten psychedelic phenethylamines in urine by gas chromatography-mass spectrometry. *J. Anal. Toxicol.* **35**(7), 459-469 (2011).

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, 05/09/2020

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