

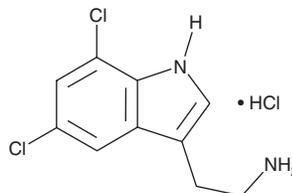
PRODUCT INFORMATION



5,7-dichloro Tryptamine (hydrochloride)

Item No. 17331

CAS Registry No.: 94850-28-9
Formal Name: 5,7-dichloro-1H-indole-3-ethanamine, monohydrochloride
MF: C₁₀H₁₀Cl₂N₂ • HCl
FW: 265.6
Purity: ≥95%
UV/Vis.: λ_{max}: 229, 294 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

5,7-dichloro Tryptamine is a ring-substituted tryptamine derivative. Although the biological effects of this compound have not been reported, tryptamines substituted with fluorine at the 5-, 6-, or 7-position have been shown to have 5-HT_{2C} receptor selectivity over both the 5-HT_{2A} and 5-HT_{2B} subtypes.^{1,2} This product is intended for forensic and research applications.

References

1. Kalir, A. and Szara, S. Synthesis and pharmacological activity of fluorinated tryptamine derivatives. *J. Med. Chem.* **6(6)**, 716-719 (1963).
2. Blair, J.B., Kurrasch-Orbaugh, D., Marona-Lewicka, D., et al. Effect of ring fluorination on the pharmacology of hallucinogenic tryptamines. *J. Med. Chem.* **43(24)**, 4701-4710 (2000).

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, 06/16/2023

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