

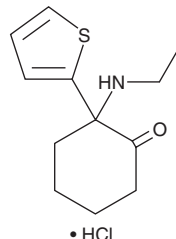
PRODUCT INFORMATION



Tiletamine (hydrochloride)

Item No. 19794

CAS Registry No.: 14176-50-2
Formal Name: 2-(ethylamino)-2-(2-thienyl)-cyclohexanone, monohydrochloride
Synonyms: CI-634, CL-399, CN 54521-2, NSC 167740
MF: C₁₂H₁₇NOS • HCl
FW: 259.8
Purity: ≥98%
UV/Vis.: λ_{max}: 236 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

Tiletamine (hydrochloride) (Item No. 19794) is an analytical reference standard that is structurally classified as an arylcyclohexylamine. It is a non-competitive NMDA receptor antagonist used in veterinary medicine for its anesthetic and sedative properties, which are comparable to ketamine (Item No. 11630).^{1,2} Abuse of this compound has been documented.² It is intended for forensic and research purposes only.

References

1. Klockgether, T., Turski, L., Schwarz, M., *et al.* Paradoxical convulsant action of a novel non-competitive N-methyl-D-aspartate (NMDA) antagonist, tiletamine. *Brain Res.* **461(2)**, 343-348 (1988).
2. Quail, M. T., Weimersheimer, P., Woolf, A. D., *et al.* Abuse of telazol: An animal tranquilizer. *J. Toxicol. Clin. Toxicol.* **39(4)**, 399-402 (2001).

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, 06/20/2023

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