PRODUCT INFORMATION



• HCI

4-chloro Pentedrone (hydrochloride)

Item No. 19383

CAS Registry No.: 2749282-78-6

1-(4-chlorophenyl)-2-(methylamino)-1-Formal Name:

pentanone, monohydrochloride

Synonyms: 4-CPD, 4-chloro-α-methylamino-Pentiophenone,

4-chloro-α-methylamino-Valerophenone

C₁₂H₁₆CINO • HCI MF:

262.2 FW: **Purity:** ≥98% Supplied as: A neat solid Storage: -20°C Stability: ≥7 years

Description

4-chloro Pentedrone (hydrochloride) (Item No. 19383) is an analytical reference material categorized as a cathinone. This product is intended for research and forensic applications.

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

Reference

1. Liu, C., Jia, W., Li, T., et al. Identification and analytical characterization of nine synthetic cathinone derivatives N-ethylhexedrone, 4-Cl-pentedrone, 4-Cl-α-EAPP, propylone, N-ethylnorpentylone, 6-MeO-bk-MDMA, α-PiHP, 4-Cl-α-PHP, and 4-F-α-PHP. Drug Test Anal. 9(8), 1162-1171 (2017).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the mater can be found on our website.

Copyright Cayman Chemical Company, 08/16/2022

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