

PRODUCT INFORMATION



α -Methylaminohexanophenone (hydrochloride)

Item No. 9002181

Formal Name: 2-(methylamino)-1-phenylhexan-1-one, monohydrochloride

Synonym: Hexedrone

MF: $C_{13}H_{19}NO \cdot HCl$

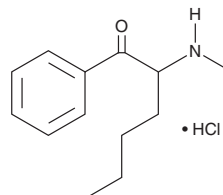
FW: 241.8

Purity: $\geq 98\%$

Supplied as: A neat solid

Storage: $-20^{\circ}C$

Stability: ≥ 2 years



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

Description

α -Methylaminohexanophenone is a substituted cathinone that is structurally similar to methcathinone (Item No. 11709), a psychoactive stimulant that is a controlled substance in the United States and other countries. It differs from methcathinone by having a butyl, rather than methyl, side chain. The physiological and toxicological properties of α -methylaminohexanophenone have not been established. This compound is intended for forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

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, 07/11/2019

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