

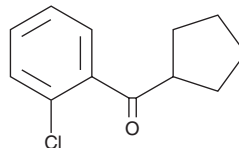
PRODUCT INFORMATION



2-Chlorophenyl Cyclopentyl Ketone

Item No. 39809

CAS Registry No.: 6740-85-8
Formal Name: (2-chlorophenyl)cyclopentyl-methanone
Synonym: o-Chlorophenyl cyclopentyl ketone
MF: C₁₂H₁₃ClO
FW: 208.7
Purity: ≥95%
Supplied as: A neat liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

2-Chlorophenyl cyclopentyl ketone is supplied as a neat liquid. A stock solution may be made by dissolving the 2-chlorophenyl cyclopentyl ketone in the solvent of choice, which should be purged with an inert gas. 2-Chlorophenyl cyclopentyl ketone is soluble in methanol.

Description

2-Chlorophenyl cyclopentyl ketone is a starting material in the synthesis of ketamine and norketamine.¹

Reference

1. Hernando, M.V., Moore, J.C., Howie, R.A., *et al.* High yielding continuous-flow synthesis of norketamine. *Org. Process Res. Dev.* **26(4)**, 1145-1151 (2022).

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, 04/29/2024

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