

PRODUCT INFORMATION



1-(4-Chlorophenyl)piperazine

Item No.14481

CAS Registry No.: 38212-33-8
Formal Name: 1-(4-chlorophenyl)-piperazine
Synonyms: 4-Chlorophenylpiperazine, *p*-Chlorophenylpiperazine, *para*-Chlorophenylpiperazine, 4-CPP, *p*-CPP, *para*-CPP, NSC 126708

MF: C₁₀H₁₃ClN₂

FW: 196.7

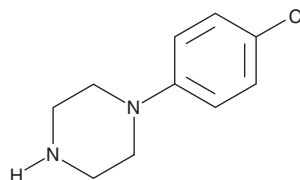
Purity: ≥98%

UV/Vis.: λ_{max}: 258 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

1-(4-Chlorophenyl)piperazine (Item No. 14481) is an analytical reference standard categorized as a piperazine. It is psychoactive and an isomer of 1-(3-chlorophenyl)piperazine (Item No. 13949).¹ This product is intended for research and forensic applications.

Reference

1. Europol-EMCDDA Joint report on a new psychoactive substance: 1-(3-chlorophenyl) piperazine (mCPP). (2005).

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