

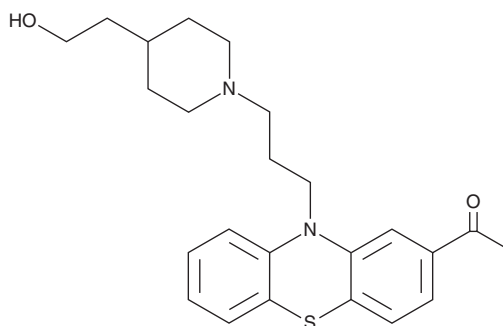
PRODUCT INFORMATION



Piperacetazine

Item No. 35220

CAS Registry No.: 3819-00-9
Formal Name: 1-[10-[3-[4-(2-hydroxyethyl)-1-piperidiny]propyl]-10H-phenothiazin-2-yl]-ethanone
Synonyms: PC-1421, SC-9794
MF: C₂₄H₃₀N₂O₂S
FW: 410.6
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 279 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Piperacetazine is supplied as a solid. A stock solution may be made by dissolving the piperacetazine in the solvent of choice, which should be purged with an inert gas. Piperacetazine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of piperacetazine in these solvents is approximately 10 and 11 mg/ml, respectively. Piperacetazine is also slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of piperacetazine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of piperacetazine in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Piperacetazine is a typical antipsychotic.¹ It inhibits spontaneous motor activity and responding for food in rats, as well as potentiates hexobarbital-induced narcosis in mice. It also inhibits lysosome acidification, as well as viral entry (IC₅₀ = <8.5 μM), in Ebola-infected Huh7 cells.² Formulations containing piperacetazine have been used in the treatment of schizophrenia.

References

1. Pitman-Moore. Piperacetazine (Quide tablets). *Clin. Pharmacol. Ther.* **10**(5), 749-754 (1969).
2. Dyall, J., Nelson, E.A., DeWald, L.E., et al. Identification of combinations of approved drugs with synergistic activity against Ebola virus in cell cultures. *J. Infect. Dis.* **218**(Suppl 5), S672-S678 (2018).

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, 12/12/2022

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