

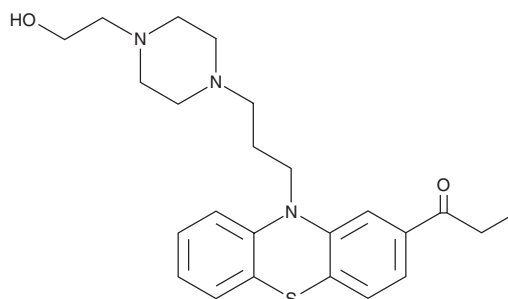
PRODUCT INFORMATION



Carfenazine

Item No. 36486

CAS Registry No.: 2622-30-2
Formal Name: 1-[10-[3-[4-(2-hydroxyethyl)-1-piperazinyl]propyl]-10H-phenothiazin-2-yl]-1-propanone
Synonyms: Carphenazine, NSC 71755
MF: C₂₄H₃₁N₃O₂S
FW: 425.6
Purity: ≥90%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Carfenazine is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Carfenazine is slightly soluble in DMSO.

Description

Carfenazine is a phenothiazine derivative.¹ It induces hypothermia in rats when administered at a dose of 10 mg/kg. Formulations containing carfenazine have previously been used in the treatment of schizophrenia.

Reference

1. Yehuda, S. and Frommer, R. The possible role of dopamine in phenothiazine-induced hypothermia in rats: An application to DA hypothesis of schizophrenia. *Int. J. Neurosci.* **7(2)**, 67-72 (1977).

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, 03/19/2026

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