

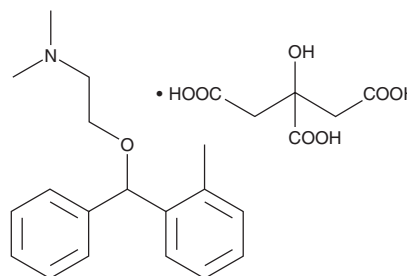
PRODUCT INFORMATION



Orphenadrine (citrate)

Item No. 39028

CAS Registry No.: 4682-36-4
Formal Name: N,N-dimethyl-2-[(2-methylphenyl)phenylmethoxy]-ethanamine, 2-hydroxy-1,2,3-propanetricarboxylate
MF: C₁₈H₂₃NO • C₆H₈O₇
FW: 461.5
Purity: ≥98%
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.

Description

Orphenadrine (citrate) (Item No. 39028) is an analytical reference standard categorized as a muscle relaxant, an anticholinergic agent, and an antihistamine.¹⁻³ Orphenadrine has been associated with misuse.^{2,3} Formulations containing orphenadrine have been used in the treatment of musculoskeletal pain. This product is intended for research and forensic applications.

References

1. Hunskaar, S. and Donnell, D. Clinical and pharmacological review of the efficacy of orphenadrine and its combination with paracetamol in painful conditions. *J. Int. Med. Res.* **19(2)**, 71-87 (1991).
2. Chiappini, S., Mosca, A., Miuli, A., et al. Misuse of anticholinergic medications: A systematic review. *Biomedicines* **10(2)**, 355 (2022).
3. Nerush, M.O., Shevyrin, V.A., Golushko, N.I., et al. Classics in chemical neuroscience: Deliriant antihistaminic drugs. *ACS Chem Neurosci.* **15(21)**, 3848-3862 (2024).

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, 12/12/2025

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