

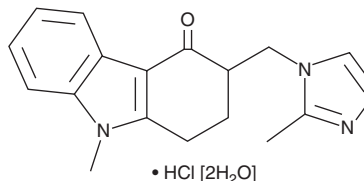
PRODUCT INFORMATION



Ondansetron (hydrochloride hydrate)

Item No. 14999

CAS Registry No.: 103639-04-9
Formal Name: 1,2,3,9-tetrahydro-9-methyl-3-[(2-methyl-1H-imidazol-1-yl)methyl]-4H-carbazol-4-one, monohydrochloride, dihydrate
MF: C₁₈H₁₉N₃O • HCl [2H₂O]
FW: 365.9
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

Ondansetron (hydrochloride hydrate) (Item No. 14999) is an analytical reference material categorized as an antiemetic.¹ Formulations containing ondansetron have been used to reduce alcohol consumption and mood disturbances in early-onset alcoholics.^{1,2} This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Hesketh, P.J. Comparative review of 5-HT₃ receptor antagonists in the treatment of acute chemotherapy-induced nausea and vomiting. *Cancer Invest.* **18(2)**, 163-173 (2000).
2. Johnson, B.A., Roache, J.D., Javors, M.A., *et al.* Ondansetron for reduction of drinking among biologically predisposed alcoholic patients: A randomized controlled trial. *JAMA* **284(8)**, 963-971 (2000).
3. Johnson, B.A., Ait-Daoud, N., Ma, J.Z., *et al.* Ondansetron reduces mood disturbance among biologically predisposed, alcohol-dependent individuals. *Alcohol Clin. Exp. Res.* **27(11)**, 1773-1779 (2003).

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, 08/26/2019

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