# **PRODUCT** INFORMATION



Methagualone

Item No. 9000636

| CAS Registry No.:  | 72-44-6  | $\sim$ |
|--|--|--------|
| Formal Name:   | 2-methyl-3-(2-methylphenyl)-4(3H)-quinazolinone  |        |
| Synonyms:  | NSC 111388, NSC 126877, NSC 631628               |        |
| MF:  | C <sub>16</sub> H <sub>14</sub> N <sub>2</sub> O |        |
| FW:  | 250.3  |        |
| Purity:  | ≥98%   |        |
| Supplied as:   | A neat solid                                     | N      |
| Storage:   | -20°C  |        |
| Stability:   | ≥1 year  |        |
| Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis. |  |        |

Description

Methaqualone (Item No. 9000636) is an analytical reference material that is structurally categorized as a quinazolinone. It is a general central nervous system depressant that has previously been approved for use as a sleep aid.<sup>1</sup> Excessive use leads to tolerance with psychological and physical dependence.<sup>2</sup> Methaqualone is regulated (Schedule I) in the United States. This product is intended for research and forensic applications.

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

## References

- 1. Ator, N. A. and Griffiths, R. R. Selectivity in the generalization profile in baboons trained to discriminate lorazepam: Benzodiazepines, barbiturates and other sedative/anxiolytics. J. Pharmacol. Exp. Ther. 282(3), 1442-1457 (1997).
- 2. Griffiths, R. R. and Johnson, M. W. Relative abuse liability of hypnotic drugs: A conceptual framework and algorithm for differentiating among compounds. J. Clin. Psychiatry 66 Suppl 9, 31-41 (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.

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, 09/30/2016

# 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