

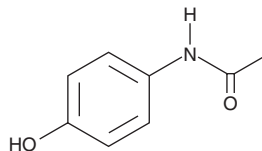
PRODUCT INFORMATION



Acetaminophen

Item No. 22629

CAS Registry No.: 103-90-2
Formal Name: N-(4-hydroxyphenyl)-acetamide
Synonyms: 4-Acetamidophenol, APAP, 4'-Hydroxyacetanilide, NSC 3991, NSC 109028, Paracetamol
MF: C₈H₉NO₂
FW: 151.2
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥6 years



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

Description

Acetaminophen (Item No. 22629) is an analytical reference material categorized as an analgesic and antipyretic agent.^{1,2} Acetaminophen has been found as an adulterant in seized heroin samples.³ Formulations containing acetaminophen have been associated with overdoses.⁴ This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 10024).

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. Vinegar, R., Truax, J.F., and Selph, J.L. Quantitative comparison of the analgesic and anti-inflammatory activities of aspirin, phenacetin and acetaminophen in rodents. *Eur. J. Pharmacol.* **37(1)**, 23-30 (1976).
2. Kobayashi, S. and Takagi, H. Fever responses to bacterial pyrogens in guinea pigs and application for screening of antipyretic agents. *Jpn. J. Pharmacol.* **18(1)**, 80-85 (1968).
3. Lurie, I.S., Driscoll, S.E., Cathapermal, S.S., et al. Determination of heroin and basic impurities for drug profiling by ultra-high-pressure liquid chromatography. *Forensic Sci. Int.* **231(1-3)**, 300-305 (2013).
4. Budnitz, D.S., Lovegrove, M.C., and Crosby, A.E. Emergency department visits for overdoses of acetaminophen-containing products. *Am. J. Prev. Med.* **40(6)**, 585-592 (2011).

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, 02/22/2021

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