

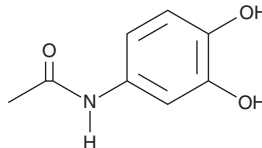
PRODUCT INFORMATION



3-Hydroxyacetaminophen

Item No. 45936

CAS Registry No.: 37519-14-5
Formal Name: N-(3,4-dihydroxyphenyl)-acetamide
Synonyms: 3-hydroxy APAP,
3'-Hydroxyacetanilide
MF: C₈H₉NO₃
FW: 167.2
Purity: ≥95%
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.

Laboratory Procedures

3-Hydroxyacetaminophen is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxyacetaminophen in the solvent of choice, which should be purged with an inert gas. 3-Hydroxyacetaminophen is soluble (≥10 mg/ml) in DMSO. 3-Hydroxyacetaminophen is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

3-Hydroxyacetaminophen is an active metabolite of acetaminophen (Item No. 10024).¹ It is formed from acetaminophen by the cytochrome P450 (CYP) isoforms CYP2C11, -2A1, -2C12, -2B1, -3A, and -1A, among others. 3-Hydroxyacetaminophen induces hepatic necrosis to a lesser extent than acetaminophen in mice.²

References

1. Harvison, P.J., Guengerich, F.P., Rashed, M.S., *et al.* Cytochrome P-450 isozyme selectivity in the oxidation of acetaminophen. *Chem. Res. Toxicol.* **1**(1), 47-52 (1998).
2. Forte, A.J., Wilson, J.M., Slattery, J.T., *et al.* The formation and toxicity of catechol metabolites of acetaminophen in mice. *Drug. Metab. Dispos.* **12**(4), 484-491 (1984).

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, 07/08/2026

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