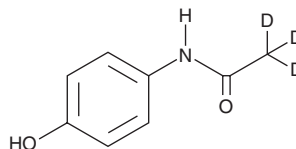


PRODUCT INFORMATION

Acetaminophen-d₃

Item No. 46351

CAS Registry No.: 60902-28-5
Formal Name: N-(4-hydroxyphenyl)-acetamide-2,2,2-d₃
Synonyms: 4-Acetamidophenol-d₃, APAP-d₃, 4'-Hydroxyacetanilide-d₃, Paracetamol-d₃
MF: C₈H₆D₃NO₂
FW: 154.2
Chemical Purity: ≥98% (Acetaminophen)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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

Acetaminophen-d₃ is intended for use as an internal standard for the quantification of acetaminophen (Item No. 10024) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Acetaminophen-d₃ is supplied as a solid. A stock solution may be made by dissolving the acetaminophen-d₃ in the solvent of choice, which should be purged with an inert gas. Acetaminophen-d₃ is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

Acetaminophen is an analgesic and antipyretic compound.^{1,2} Unlike many NSAIDs, which inhibit both COX-1 and COX-2, early studies suggested that acetaminophen is a poor inhibitor of both isoforms.^{3,4} However, it does inhibit COX-2 by 83% and COX-1 by 56% in human blood *ex vivo*, albeit at a high 1,000 mg dose, with IC₅₀ values of 25.8 and 113.7 μM, respectively.⁵ Acetaminophen is enzymatically and non-enzymatically converted to several reactive metabolites that contribute to adverse or indirect effects, including liver injury.⁶⁻⁸ At toxic doses, the acetaminophen metabolite N-acetyl-4-benzoquinone imine (NAPQI; Item No. 16115) depletes glutathione reserves in the liver, leading to an accumulation of NAPQI and subsequent hepatocyte necrosis.⁹ Acetaminophen decreases glutathione levels and reduces glutathione peroxidase activity in mice when administered at a dose of 250 mg/kg and induces ferroptotic cell death in primary mouse hepatocytes, an effect that can be blocked by the ferroptosis inhibitor ferrostatin-1 (Item No. 17729).^{10,11} Acetaminophen has analgesic and antipyretic properties in animal models.^{1,2}

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. Mitchell, J.A., Akarasereenont, P., Thiemermann, C., *et al.* Selectivity of nonsteroidal antiinflammatory drugs as inhibitors of constitutive and inducible cyclooxygenase. *Proc. Natl. Acad. Sci. USA* **90(24)**, 11693-11697 (1993).

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, 08/28/2026

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

PRODUCT INFORMATION

4. Warner, T.D., Giuliano, F., Vojnovic, I., *et al.* Nonsteroid drug selectivities for cyclo-oxygenase-1 rather than cyclo-oxygenase-2 are associated with human gastrointestinal toxicity: A full *in vitro* analysis. *Proc. Natl. Acad. Sci. USA* **96**(13), 7563-7568 (1999).
5. Hinz, B., Cheremina, O., and Brune, K. Acetaminophen (paracetamol) is a selective cyclooxygenase-2-inhibitor in man. *FASEB J.* **22**(2), 383-390 (2007).
6. Hinson, J.A. Reactive metabolites of phenacetin and acetaminophen: A review. *Environ. Health Perspect.* **49**, 71-79 (1983).
7. Högestätt, E.D., Jönsson, B.A.G., Ermund, A., *et al.* Conversion of acetaminophen to the biactive N-acylphenolamine AM404 via fatty acid amide hydrolase-dependent arachidonic acid conjugation in the nervous system. *J. Biol. Chem.* **280**(36), 31405-31412 (2005).
8. Han, D., Dara, L., Win, S., *et al.* Regulation of drug-induced liver injury by signal transduction pathways: Critical role of mitochondria. *Trends Pharmacol. Sci.* **34**(4), 243-253 (2013).
9. Qiu, Y., Benet, L.Z., and Burlingame, A.L. Identification of the hepatic protein targets of reactive metabolites of acetaminophen *in vivo* in mice using two-dimensional gel electrophoresis and mass spectrometry. *J. Biol. Chem.* **273**(28), 17940-17953 (1998).
10. Tirmenstein, M.A. and Nelson, S.D. Acetaminophen-induced oxidation of protein thiols. Contribution of impaired thiol-metabolizing enzymes and the breakdown of adenine nucleotides. *The Journal of Biological Chemistry* **265**(6), 3059-3065 (1990).
11. Lőrincz, T., Jemnitz, K., Kardon, T., *et al.* Ferroptosis is involved in acetaminophen induced cell death. *Pathol. Oncol. Res.* **21**(4), 1115-1121 (2015).

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