

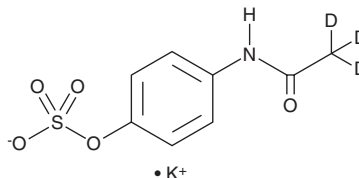
PRODUCT INFORMATION



Acetaminophen-d₃ sulfate (potassium salt)

Item No. 46350

CAS Registry No.: 1188263-45-7
Formal Name: N-[4-(sulfooxy)phenyl]-acetamide-2,2,2-d₃,
monopotassium salt
Synonyms: 4-Acetamidophenol-d₃ sulfate,
APAP-d₃ sulfate,
4'-Hydroxyacetanilide-d₃ sulfate,
Paracetamol-d₃ sulfate
MF: C₈H₅D₃NO₅S • K
FW: 272.3
Chemical Purity: ≥98% (Acetaminophen sulfate (potassium salt))
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 sulfate (potassium salt) is intended for use as an internal standard for the quantification of acetaminophen sulfate (Item No. 22571) 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 sulfate (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the acetaminophen-d₃ sulfate (potassium salt) in the solvent of choice, which should be purged with an inert gas. Acetaminophen-d₃ sulfate (potassium salt) is sparingly soluble (1-10 mg/ml) in acetonitrile.

Description

Acetaminophen sulfate is an inactive metabolite of the antipyretic and analgesic agent acetaminophen (Item No. 10024).¹ It is formed from acetaminophen via the sulfotransferase (SULT) isoforms SULT2A1, SULT1E1, SULT1A1, SULT1A3, and SULT1A4.

Reference

1. Mazaleuskaya, L.L., Sangkuhl, K., Thorn, C.F., *et al.* PharmGKB summary: Pathways of acetaminophen metabolism at the therapeutic versus toxic doses. *Pharmacogenet. Genomics* **25**(8), 416-426 (2015).

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, 09/21/2026

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