

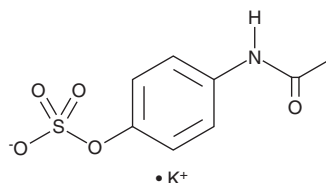
# PRODUCT INFORMATION



## Acetaminophen sulfate (potassium salt)

Item No. 22571

**CAS Registry No.:** 32113-41-0  
**Formal Name:** N-[4-(sulfooxy)phenyl]-acetamide, monopotassium salt  
**Synonyms:** 4-Acetamidophenol sulfate, APAP sulfate, 4'-Hydroxyacetanilide sulfate, Paracetamol sulfate  
**MF:** C<sub>8</sub>H<sub>8</sub>NO<sub>5</sub>S • K  
**FW:** 269.3  
**Purity:** ≥98%  
**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 sulfate (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the acetaminophen sulfate (potassium salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

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

### References

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, 11/21/2022

#### 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