

# PRODUCT INFORMATION



## Acetaminophen Glucuronide (sodium salt)

Item No. 27850

CAS Registry No.: 120595-80-4

Formal Name: 4-(acetylamino)phenyl β-D-glucopyranosiduronic acid, monosodium salt

Synonyms: 4-Acetamidophenol Glucuronide, APAP Glucuronide, 4'-Hydroxyacetanilide Glucuronide, Paracetamol Glucuronide

MF: C<sub>14</sub>H<sub>16</sub>NO<sub>8</sub> • Na

FW: 349.3

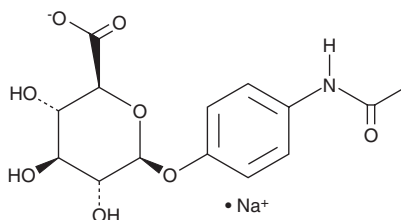
Purity: ≥98%

UV/Vis.: λ<sub>max</sub>: 246 nm

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 glucuronide (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the acetaminophen glucuronide (sodium salt) in the solvent of choice, which should be purged with an inert gas. Acetaminophen glucuronide (sodium salt) is slightly soluble in methanol.

Acetaminophen glucuronide (sodium salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Acetaminophen glucuronide is an inactive metabolite of the analgesic and antipyretic agent acetaminophen (Item No. 10024).<sup>1</sup> It is formed *via* glucuronidation of acetaminophen by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A6, UGT1A9, UGT1A1, and UGT2B15.

### 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, 11/14/2022

#### 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