

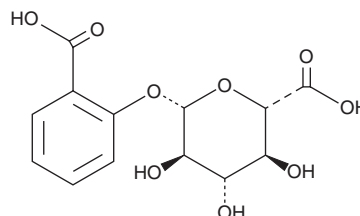
# PRODUCT INFORMATION



## 1-Salicylate Glucuronide

Item No. 34820

**CAS Registry No.:** 7695-70-7  
**Formal Name:** 2-carboxyphenyl β-D-glucopyranosiduronic acid  
**Synonyms:** Salicyl Phenolic Glucuronide, Salicylic Acid Phenolic Glucuronide  
**MF:** C<sub>13</sub>H<sub>14</sub>O<sub>9</sub>  
**FW:** 314.2  
**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

1-Salicylate glucuronide is supplied as a solid. A stock solution may be made by dissolving the 1-salicylate glucuronide in the solvent of choice, which should be purged with an inert gas. 1-Salicylate glucuronide is slightly soluble in methanol.

1-Salicylate glucuronide 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

1-Salicylate glucuronide is a metabolite of salicylic acid and aspirin (Item No. 70260).<sup>1</sup> It is formed from salicylic acid primarily by the UDP-glucuronosyltransferase (UGT) isoform UGT1A9 but also by a variety of other UGT isoforms and from aspirin *via* salicylic acid as an intermediate.

### Reference

1. Kuehl, G.E., Bigler, J., Potter, J.D., *et al.* Glucuronidation of the aspirin metabolite salicylic acid by expressed UDP-glucuronosyltransferases and human liver microsomes. *Drug Metab. Dispos.* **34**(2), 199-202 (2006).

#### 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, 12/13/2022

#### 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