

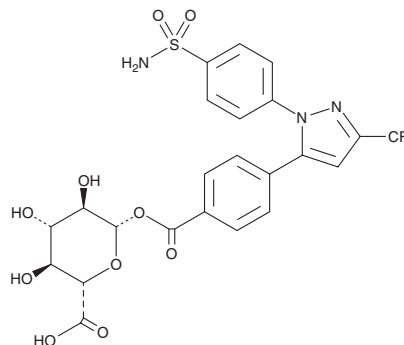
# PRODUCT INFORMATION



## Celecoxib Carboxylic Acid Acyl-β-D-Glucuronide

Item No. 22457

**CAS Registry No.:** 264236-79-5  
**Formal Name:** β-D-glucopyranuronic acid, 1-[4-[1-[4-(aminosulfonyl)phenyl]-3-(trifluoromethyl)-1H-pyrazol-5-yl]benzoate]  
**MF:** C<sub>23</sub>H<sub>20</sub>F<sub>3</sub>N<sub>3</sub>O<sub>10</sub>S  
**FW:** 587.5  
**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

Celecoxib carboxylic acid acyl-β-D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the celecoxib carboxylic acid acyl-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Celecoxib carboxylic acid acyl-β-D-glucuronide is soluble in DMSO. Celecoxib carboxylic acid acyl-β-D-glucuronide is slightly soluble in methanol.

### Description

Celecoxib carboxylic acid acyl-β-D-glucuronide is a metabolite of the non-steroidal anti-inflammatory drug (NSAID) and COX-2 inhibitor celecoxib (Item No. 10008672).<sup>1</sup> It is formed *via* a multi-step process in which celecoxib undergoes hydroxylation to form hydroxy celecoxib (Item No. 33994) by the cytochrome P450 (CYP) isoforms CYP2C9, CYP3A4, and CYP2D6, followed by oxidation to form celecoxib carboxylic acid (Item No. 34062), and then conjugation to the β-glucuronide.<sup>2,3</sup>

### References

1. Paulson, S.K., Hribar, J.D., Liu, N.W., *et al.* Metabolism and excretion of [<sup>14</sup>C]celecoxib in healthy male volunteers. *Drug Metab. Dispos.* **28(3)**, 308-314 (2000).
2. Ma, Y., Gao, S., and Hu, M. Quantitation of celecoxib and four of its metabolites in rat blood by UPLC-MS/MS clarifies their blood distribution patterns and provides more accurate pharmacokinetics profiles. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **1001**, 202-211 (2015).
3. Siu, Y.A., Hao, M.-H., Dixit, V., *et al.* Celecoxib is a substrate of CYP2D6: Impact on celecoxib metabolism in individuals with CYP2C9\*3 variants. *Drug Metab. Pharmacokinet.* **33(5)**, 219-227 (2018).

#### 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, 10/17/2022

#### 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