

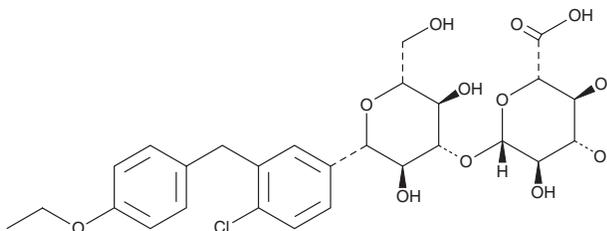
# PRODUCT INFORMATION



## Dapagliflozin-3-O-β-D-Glucuronide

Item No. 22458

**CAS Registry No.:** 1351438-75-9  
**Formal Name:** (1S)-1,5-anhydro-1-C-[4-chloro-3-[(4-ethoxyphenyl)methyl]phenyl]-3-O-β-D-glucopyranuronosyl-D-glucitol  
**MF:** C<sub>27</sub>H<sub>33</sub>ClO<sub>12</sub>  
**FW:** 585.0  
**Purity:** ≥95%  
**Supplied as:** A crystalline 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

Dapagliflozin-3-O-β-D-glucuronide is supplied as a crystalline solid. A stock solution may be made by dissolving the dapagliflozin-3-O-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Dapagliflozin-3-O-β-D-glucuronide is soluble in organic solvents such as methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Dapagliflozin-3-O-β-D-glucuronide is a metabolite of dapagliflozin (Item No. 11574). Dapagliflozin is a first generation selective sodium glucose cotransporter (SGLT) 2 inhibitor that blocks glucose transport.<sup>1</sup> Dapagliflozin is metabolized by uridine diphosphate glucuronosyltransferase (UGT) 1A9 to dapagliflozin-3-O-β-D-glucuronide, which is 2,600-fold less potent than the parent compound with regard to SGLT2 inhibition.<sup>2</sup>

### References

1. Chao, E.C. and Henry, R.R. SGLT2 inhibition - A novel strategy for diabetes treatment. *Nat. Rev. Drug Discov.* **9**(7), 551-559 (2010).
2. van der Walt, J.-S., Hong, Y., Zhang, L., *et al.* A nonlinear mixed effects pharmacokinetic model for dapagliflozin and dapagliflozin 3-O-glucuronide in renal or hepatic impairment. *CPT Pharmacometrics Syst. Pharmacol.* **2**(e42), 1-9 (2013).

#### 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, 11/07/2022

#### 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