

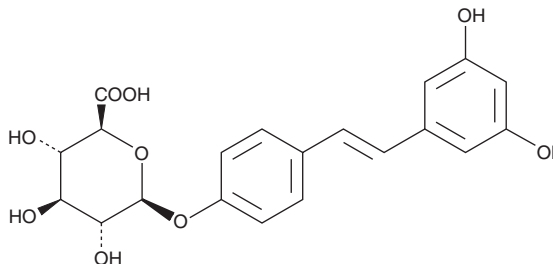
# PRODUCT INFORMATION



## *trans*-Resveratrol-4'-O-D-Glucuronide

Item No. 13833

**CAS Registry No.:** 387372-20-5  
**Formal Name:** 4-[(1E)-2-(3,5-dihydroxyphenyl)ethenyl]phenyl-β-D-glucopyranosiduronic acid  
**Synonym:** *trans*-R4'G  
**MF:** C<sub>20</sub>H<sub>20</sub>O<sub>9</sub>  
**FW:** 404.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 213, 303 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

*trans*-Resveratrol-4'-O-D-glucuronide is supplied as a crystalline solid. A stock solution may be made by dissolving the *trans*-resveratrol-4'-O-D-glucuronide in the solvent of choice. *trans*-Resveratrol-4'-O-D-glucuronide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of *trans*-resveratrol-4'-O-D-glucuronide in these solvents is approximately 0.2, 10, and 15 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of *trans*-resveratrol-4'-O-D-glucuronide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of *trans*-resveratrol-4'-O-D-glucuronide in PBS, pH 7.2, is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

*trans*-Resveratrol-4'-O-D-glucuronide is a phase II metabolite of the antioxidant *trans*-resveratrol (Item No. 70675).<sup>1</sup> *trans*-Resveratrol-4'-O-D-glucuronide is formed *via* glucuronidation of *trans*-resveratrol by the UDP-glucuronosyltransferase (UGT) isoform UGT1A9.

### Reference

1. Matin, B., Sherbini, A.A., Alam, N., *et al.* Resveratrol glucuronidation *in vitro*: Potential implications of inhibition by probenecid. *J. Pharm. Pharmacol.* (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, 01/09/2019

### 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