

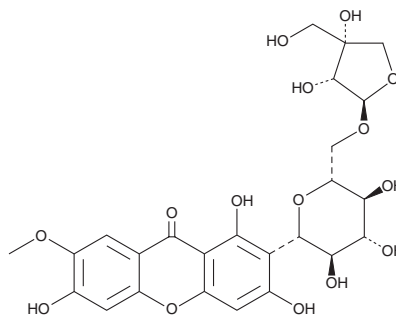
# PRODUCT INFORMATION



## Polygalaxanthone III

Item No. 30621

**CAS Registry No.:** 162857-78-5  
**Formal Name:** 2-(6-O-D-apio-β-D-furanosyl-β-D-glucopyranosyl)-1,3,6-trihydroxy-7-methoxy-9H-xanthen-9-one  
**MF:** C<sub>25</sub>H<sub>28</sub>O<sub>15</sub>  
**FW:** 568.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 242, 258, 316, 360 nm  
**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

Polygalaxanthone III is supplied as a crystalline solid. A stock solution may be made by dissolving the polygalaxanthone III in the solvent of choice, which should be purged with an inert gas. Polygalaxanthone III is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of polygalaxanthone III in these solvents is approximately 15 and 10 mg/ml, respectively.

Polygalaxanthone III is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, polygalaxanthone III should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Polygalaxanthone III has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Polygalaxanthone III is a xanthone glycoside that has been found in *P. hongkongensis* and has antioxidant activity.<sup>1</sup> It scavenges hydroxyl and 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in cell-free assays (IC<sub>50</sub>s = 83.5 and 76.1 μM, respectively).

### Reference

1. Wu, J.-F., Chen, S.-B., Gao, J.-C., *et al.* Xanthone glycosides from herbs of *Polygala hongkongensis* Hemsl and their antioxidant activities. *J. Asian Nat. Prod. Res.* **10(7-8)**, 673-678 (2008).

#### 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, 12/12/2022

#### 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