

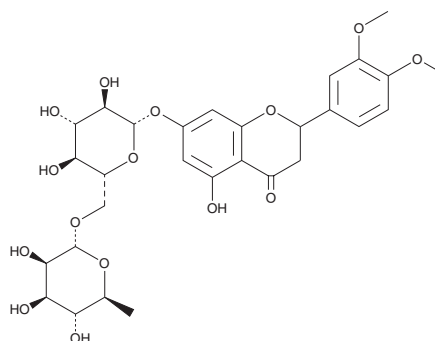
# PRODUCT INFORMATION



## Methylhesperidin

Item No. 30049

**CAS Registry No.:** 1616116-58-5  
**Formal Name:** 7-[[6-O-(6-deoxy- $\alpha$ -L-mannopyranosyl)- $\beta$ -D-glucopyranosyl]oxy]-2-(3,4-dimethoxyphenyl)-2,3-dihydro-5-hydroxy-4H-1-benzopyran-4-one  
**MF:** C<sub>29</sub>H<sub>36</sub>O<sub>15</sub>  
**FW:** 624.6  
**Purity:**  $\geq 90\%$  (mixture of isomers)  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 226, 282 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Methylhesperidin is supplied as a crystalline solid. A stock solution may be made by dissolving the methylhesperidin in the solvent of choice, which should be purged with an inert gas. Methylhesperidin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of methylhesperidin in these solvents is approximately 5 and 3 mg/ml, respectively.

Methylhesperidin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, methylhesperidin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Methylhesperidin has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Methylhesperidin is a methylated form of the flavanone rutoside hesperidin (Item No. 18646).<sup>1</sup>

### Reference

1. Li, J. and Wang, S. Molecular spectroscopic on interaction between Methyl hesperidin and Buman serum albumin. *Spectrochim. Acta A Mol. Biomol. Spectrosc.* **102**, 200-204 (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, 11/14/2022

#### 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](http://WWW.CAYMANCHEM.COM)