

# PRODUCT INFORMATION



## Methylophiopogonanone A

Item No. 30192

CAS Registry No.: 74805-92-8  
Formal Name: (3R)-3-(1,3-benzodioxol-5-ylmethyl)-2,3-dihydro-5,7-dihydroxy-6,8-dimethyl-4H-1-benzopyran-4-one

MF: C<sub>19</sub>H<sub>18</sub>O<sub>6</sub>

FW: 342.3

Purity: ≥98%

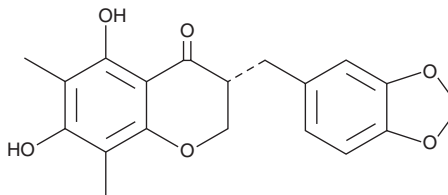
UV/Vis.: λ<sub>max</sub>: 230 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/Ophiopogonis radix



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

### Laboratory Procedures

Methylophiopogonanone A is supplied as a crystalline solid. A stock solution may be made by dissolving the methylophiopogonanone A in the solvent of choice, which should be purged with an inert gas. Methylophiopogonanone A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of methylophiopogonanone A in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Methylophiopogonanone A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, methylophiopogonanone A should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Methylophiopogonanone A has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Methylophiopogonanone A is a homoisoflavonoid that has been found in *O. japonicus* roots and has diverse biological activities.<sup>1-4</sup> It increases the uptake of rosuvastatin (Item No. 18813) and atorvastatin (Item No. 10493) in HEK293T cells expressing human organic anion transporting polypeptide 1B1 (OATP1B1; EC<sub>50</sub>s = 11.33 and 6 μM, respectively).<sup>2</sup> Methylophiopogonanone A scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) and ABTS (Item No. 27317) radicals in cell-free assays.<sup>1</sup> *In vivo*, methylophiopogonanone A (10 mg/kg) decreases body weight gain and reduces serum and hepatic lipid levels in a rat model of high-fat diet-induced hyperlipidemia.<sup>3</sup> It also reduces infarct size and myocardial apoptosis in mouse model of ischemia-reperfusion injury induced by transient occlusion of the left anterior descending coronary artery.<sup>4</sup>

### References

1. Wang, Y., Liu, F., Liang, Z., et al. *Iran J. Pharm. Res.* **16**(1), 357-365 (2017).
2. Chen, L., Liu, L., Chen, Y., et al. *Xenobiotica* **49**(10), 1221-1228 (2019).
3. Li, Z., Wu, Y.Y., and Yu, B.X. *Braz. J. Med. Biol. Res.* **53**(3), e9201 (2020).
4. He, F., Xu, B.L., Chen, C., et al. *Acta. Pharmacol. Sin.* **37**(6), 763-771 (2016).

#### 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, 10/20/2025

#### 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