

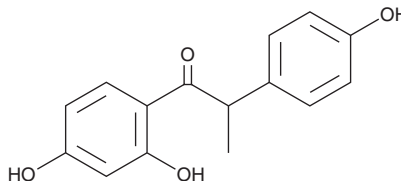
PRODUCT INFORMATION



O-Desmethylangolensin

Item No. 36435

CAS Registry No.: 21255-69-6
Formal Name: 1-(2,4-dihydroxyphenyl)-2-(4-hydroxyphenyl)-1-propanone
Synonym: O-DMA
MF: C₁₅H₁₄O₄
FW: 258.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

O-Desmethylangolensin (O-DMA) is supplied as a solid. A stock solution may be made by dissolving the O-DMA in the solvent of choice, which should be purged with an inert gas. O-DMA is soluble in DMSO and methanol.

Description

O-DMA is a phytoestrogen and an active metabolite of daidzein (Item No. 10005166).¹ It is formed from daidzein by gut microbiota. O-DMA (2.5-50 μM) is cytotoxic to MDA-MB-231 cells.²

References

1. Frankenfeld, C.L. O-desmethylangolensin: The importance of equol's lesser known cousin to human health. *Adv. Nutr.* **2(4)**, 317-324 (2011).
2. Magee, P.J., McGlynn, H., and Rowland, I.R. Differential effects of isoflavones and lignans on invasiveness of MDA-MB-231 breast cancer cells in vitro. *Cancer Lett.* **208(1)**, 35-41 (2004).

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, 10/25/2022

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