

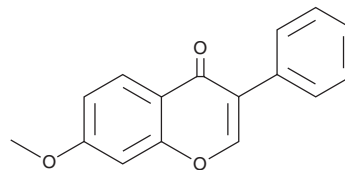
# PRODUCT INFORMATION



## 7-Methoxyisoflavone

Item No. 33419

**CAS Registry No.:** 1621-56-3  
**Formal Name:** 7-methoxy-3-phenyl-4H-1-benzopyran-4-one  
**Synonym:** 5-methyl-7-Methoxyisoflavone  
**MF:** C<sub>16</sub>H<sub>12</sub>O<sub>3</sub>  
**FW:** 252.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250 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

7-Methoxyisoflavone is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-methoxyisoflavone in the solvent of choice, which should be purged with an inert gas. 7-Methoxyisoflavone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 7-methoxyisoflavone in these solvents is approximately 10 and 30 mg/ml, respectively.

### Description

7-Methoxyisoflavone is a synthetic isoflavone and an isomer of 7-methoxyflavone (Item No. 33497).<sup>1,2</sup> It inhibits mitosis induced by concanavalin A (Item No. 14951) in human peripheral blood mononuclear cells (PBMCs) with an IC<sub>50</sub> value of 1.36 μg/ml.<sup>1</sup> 7-Methoxyisoflavone (10, 20, and 40 μM) inhibits the growth of HCT116 colon cancer cells in a concentration-dependent manner.<sup>2</sup>

### References

1. Hirano, T., Oka, K., Kawashima, E., *et al.* Effects of synthetic and naturally occurring flavonoids on mitogen-induced proliferation of human peripheral-blood lymphocytes. *Life Sci.* **45(15)**, 1407-1411 (1989).
2. Hyun, J., Shin, S.Y., So, K.M., *et al.* Isoflavones inhibit the clonogenicity of human colon cancer cells. *Bioorg. Med. Chem. Lett.* **22(8)**, 2664-2669 (2012).

#### 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, 11/07/2022

#### 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