

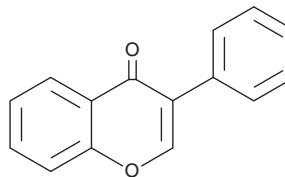
PRODUCT INFORMATION



Isoflavone

Item No. 33964

CAS Registry No.: 574-12-9
Formal Name: 3-phenyl-4H-1-benzopyran-4-one
Synonyms: 3-Phenylchromone, NSC 135405
MF: $C_{15}H_{10}O_2$
FW: 222.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Isoflavone is supplied as a solid. A stock solution may be made by dissolving the isoflavone in the solvent of choice, which should be purged with an inert gas. Isoflavone is slightly soluble in chloroform and methanol.

Description

Isoflavone is a polyketide synthase-derived flavonoid that has been found in *G. max*.^{1,2} It is the scaffold of several phytoestrogens including genistein (Item No. 10005167), daidzein (Item No. 10005166), and glycitein (Item No. 14162).

References

1. Austin, M.B. and Noel, J.P. The chalcone synthase superfamily of type III polyketide synthases. *Nat. Prod. Rep.* **20**(1), 79-110 (2003).
2. Křížová, L., Dadáková, K., Kašparovská, J., *et al.* Isoflavones. *Molecules* **24**(6), 1076 (2019).

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