

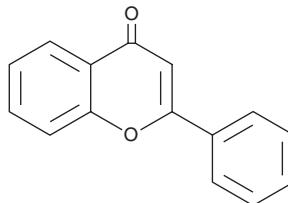
PRODUCT INFORMATION



Flavone

Item No. 34028

CAS Registry No.: 525-82-6
Formal Name: 2-phenyl-4H-1-benzopyran-4-one
Synonyms: NSC 19028, 2-Phenylchormone
MF: C₁₅H₁₀O₂
FW: 222.2
Purity: ≥98%
UV/Vis.: λ_{max}: 251, 296 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/Unidentified sp.



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

Laboratory Procedures

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

Flavone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, flavone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Flavone has a solubility of approximately 0.14 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Flavone is a polyketide synthase-derived polyphenolic flavonoid that has been found in various edible plants.¹ It is the backbone scaffold of many pharmacologically active flavonoids, including luteolin (Item No. 10004161), apigenin (Item No. 10010275), and wogonin (Item No. 14248).^{1,2} Flavone reduces proliferation of (EC₅₀ = 54.8 μM), and induces apoptosis in, HT-29 colon cancer cells.¹ It enhances dengue virus type 2 (DENV-2) infectivity in Vero cells when used at a concentration of 100 μg/ml.²

References

1. Wenzel, U., Kuntz, S., Brendel, M.D., *et al.* Dietary flavone is a potent apoptosis inducer in human colon carcinoma cells. *Cancer Res.* **60**(14), 3823-3831 (2000).
2. Zandi, K., Lani, R., Wong, P.-F., *et al.* Flavone enhances dengue virus type-2 (NGC strain) infectivity and replication in Vero cells. *Molecules* **17**(3), 2437-2445 (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, 12/08/2022

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