

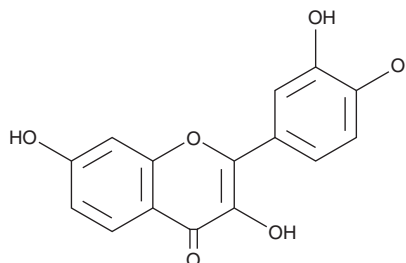
PRODUCT INFORMATION



Fisetin

Item No. 15246

CAS Registry No.: 528-48-3
Formal Name: 2-(3,4-dihydroxyphenyl)-3,7-dihydroxy-4H-1-benzopyran-4-one
Synonyms: CI-75620, NSC 407010, NSC 656275
MF: C₁₅H₁₀O₆
FW: 286.2
Purity: ≥90%
UV/Vis.: λ_{max}: 250, 320, 364 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

Fisetin is supplied as a crystalline solid. A stock solution may be made by dissolving the fisetin in the solvent of choice, which should be purged with an inert gas. Fisetin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fisetin in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Fisetin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fisetin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fisetin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fisetin is a flavonoid that has been found in various fruits and vegetables and has diverse biological activities.¹⁻⁴ It scavenges radicals in a Trolox equivalent antioxidant capacity (TEAC) assay and inhibits oxidative burst induced by N-formyl-methionyl-leucyl-phenylalanine (fMLP; Item No. 21495) in isolated human polymorphonuclear neutrophils (PMNs) when used at a concentration of 10 μM.^{1,3} Fisetin (10-60 μM) induces cell cycle arrest at the G₁ phase and apoptosis in, as well as inhibits the proliferation of, LNCaP prostate cancer cells.² It inhibits LPS-induced production of nitrite, prostaglandin E₂ (PGE₂; Item No. 14010), and NF-κB activation in RAW 264.7 cells.³ Fisetin (1 mg/animal) reduces tumor growth in a CWR22Ru1 prostate cancer mouse xenograft model.⁴

References

1. Khan, N., Syed, D.N., Ahmad, N., *et al.* Fisetin: A dietary antioxidant for health promotion. *Antioxid. Redox Signal.* **19(2)**, 151-162 (2013).
2. Khan, N., Afaq, F., Syed, D.N., *et al.* Fisetin, a novel dietary flavonoid, causes apoptosis and cell cycle arrest in human prostate cancer LNCaP cells. *Carcinogenesis* **29(5)**, 1049-1056 (2008).
3. Wang, L., Tu, Y.-C., Lian, T.-W., *et al.* Distinctive antioxidant and antiinflammatory effects of flavonols. *J. Agric. Food Chem.* **54(26)**, 9798-9804 (2006).
4. Khan, N., Asim, M., Afaq, F., *et al.* A novel dietary flavonoid fisetin inhibits androgen receptor signaling and tumor growth in athymic nude mice. *Cancer Res.* **68(20)**, 8555-8563 (2008).

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, 12/06/2022

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