

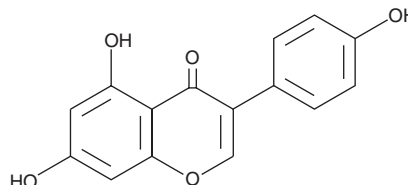
PRODUCT INFORMATION



Genistein

Item No. 10005167

CAS Registry No.: 446-72-0
Formal Name: 5,7-dihydroxy-3-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
Synonyms: CI-75610, NSC 36586
MF: C₁₅H₁₀O₅
FW: 270.2
Purity: ≥98%
UV/Vis.: λ_{max}: 261 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Styphnolobium japonicum*



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

Laboratory Procedures

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

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

Description

Genistein is an isoflavonoid phytoestrogen that has been found in soybeans (*G. max/S. hispida*) and has kinase inhibitory, anticancer, pro-cancer, hepatoprotective, and antiviral properties.¹ It inhibits the tyrosine kinases EGFR, pp50^{v-Src}, and pp110^{gag-fes} (IC₅₀s = 6, 7-8, and 6.5 μg/ml, respectively) and decreases EGF-induced serine, threonine, and tyrosine phosphorylation of EGFR in A431 cells when used at a concentration of 20 μg/ml.² Genistein inhibits proliferation and induces apoptosis in a variety of cancer cells, including Bel 7402 hepatocellular carcinoma cells when used at a concentration of 10 μg/ml.^{1,3} It reduces tumor invasion and angiogenesis in a Bel 7402 mouse subrenal capsule xenograft model when administered at a dose of 50 mg/kg per day.³ However, when administered at the same dose on postnatal days 1-5, genistein increases the incidence of uterine adenocarcinoma in a mouse model of cancer induced by the estrogen receptor agonist diethylstilbestrol (DES; Item No. 10006876).⁴ It reduces lipid accumulation and inflammation in the liver of ovariectomized (OVX) and non-OVX female rats in a model of high-fat high-fructose diet-induced nonalcoholic hepatosteatosis (NASH) when administered at a dose of 16 mg/kg per day.⁵ Genistein (10 μM) also inhibits HIV-1 DNA synthesis in resting CD4⁺ T cells.⁶

References

1. Spagnuolo, C., Russo, G.L., Orhan, I.E., et al. *Adv. Nutr.* **6(4)**, 408-419 (2015).
2. Akiyama, T., Ishida, J., Nakagawa, S., et al. *J. Biol. Chem.* **262(1)**, 5592-5595 (1987).
3. Gu, Y., Zhu, C.-F., Iwamoto, H., et al. *World J. Gastroenterol.* **11(41)**, 6512-6517 (2005).
4. Newbold, R.R., Banks, E.P., Bullock, B., et al. *Cancer Res.* **61**, 4325-4328 (2001).
5. Pummoung, S., Werawatganon, D., Klaikeaw, N., et al. *Pharmacogn. Mag.* **14(55)**, 20-24 (2018).
6. Guo, J., Xu, X., Rasheed, T.K., et al. Genistein interferes with SDF-1- and HIV-mediated actin dynamics and inhibits HIV infection of resting CD4 T cells. *Retrovirology* **10**, 62 (2013).

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, 11/04/2022

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