

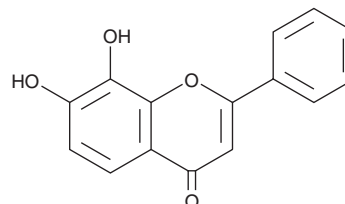
PRODUCT INFORMATION



7,8-Dihydroxyflavone

Item No. 16946

CAS Registry No.: 38183-03-8
Formal Name: 7,8-dihydroxy-2-phenyl-4H-1-benzopyran-4-one
Synonym: 7,8-DHF
MF: C₁₅H₁₀O₄
FW: 254.2
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 270 nm
Supplied as: A crystalline solid
Storage: -20°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

7,8-Dihydroxyflavone (7,8-DHF) is supplied as a crystalline solid. A stock solution may be made by dissolving the 7,8-DHF in the solvent of choice, which should be purged with an inert gas. 7,8-DHF is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 7,8-DHF in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

7,8-DHF is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 7,8-DHF should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 7,8-DHF has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

7,8-DHF is a flavone that has been found in *G. aesculifolia* and has diverse biological activities.¹⁻⁴ It binds to tropomyosin-related kinase B (TrkB; K_d = 320 nM) and induces TrkB autophosphorylation in primary mouse hippocampal neurons.¹ *In vivo*, 7,8-DHF (5 mg/kg) reduces neuronal apoptosis and reduces infarct volume in a mouse model of ischemia-reperfusion injury induced by middle cerebral artery occlusion (MCAO). It reduces striatal nigral neuronal apoptosis in a mouse model of MPTP-induced Parkinson's disease. 7,8-DHF (5 mg/kg) increases survival, neocortical volume, and whole brain volume in an N171-82Q mouse model of Huntington's disease.² It increases bone mineral density and reduces bone loss in a rat model of ovariectomy-induced osteoporosis when administered at a dose of 10 mg/kg.³ 7,8-DHF (5 mg/kg per day, i.p.) reduces blood glucose levels, pancreatic vascular congestion, and inflammatory cell infiltration, as well as hepatic vascular congestion and DNA damage-inducible transcript 3 (DDIT3) and glucose-regulated protein 78 (GRP78) levels, markers of endoplasmic reticulum (ER) stress, in a mouse model of cafeteria diet-induced metabolic syndrome.⁴

References

1. Jang, S.W., Liu, X., Yepes, M., *et al.* A selective TrkB agonist with potent neurotrophic activities by 7,8-dihydroxyflavone. *Proc. Natl. Acad. Sci. USA* **107**(6), 2687-2692 (2010).
2. Schindler, U., Strobel, H., Schönafinger, K., *et al.* Biochemistry and pharmacology of novel anthranilic acid derivatives activating heme-oxidized soluble guanylyl cyclase. *Mol. Pharmacol.* **69**(4), 1260-1268 (2006).
3. Xue, F., Zhao, Z., Gu, Y., *et al.* 7,8-Dihydroxyflavone modulates bone formation and resorption and ameliorates ovariectomy-induced osteoporosis. *Elife* **10**, 64872 (2021).
4. Sahin, E., Saglam, N., Erdem, S., *et al.* 7,8-Dihydroxyflavone alleviates endoplasmic reticulum stress in cafeteria diet-induced metabolic syndrome. *Life Sci.* **306**, 120781 (2022).

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, 10/17/2023

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