

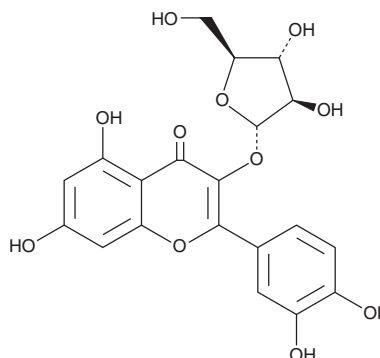
PRODUCT INFORMATION



Avicularin

Item No. 33918

CAS Registry No.: 572-30-5
Formal Name: 3-(α -L-arabinofuranosyloxy)-2-(3,4-dihydroxyphenyl)-5,7-dihydroxy-4H-1-benzopyran-4-one
Synonym: Quercetin 3-O- α -L-arabinofuranoside
MF: C₂₀H₁₈O₁₁
FW: 434.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 258, 359 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Avicularin is supplied as a solid. A stock solution may be made by dissolving the avicularin in the solvent of choice, which should be purged with an inert gas. Avicularin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of avicularin in these solvents is approximately 5, 30, and 50 mg/ml, respectively.

Description

Avicularin is a flavonoid glycoside that has been found in *R. sachalinensis* and has diverse biological activities.¹⁻⁵ It inhibits fatty acid synthase (FASN) and aldose reductase (IC₅₀s = 6.15 and 19.05 μ M for the chicken liver and human enzymes, respectively).^{1,2} Avicularin scavenges DPPH and superoxide radicals, as well as inhibits copper-induced oxidation of LDL in cell-free assays (IC₅₀s = 64.3, 6, and 3.8 μ M, respectively).³ It inhibits LPS-induced increases in nitric oxide (NO) and prostaglandin E₂ (PGE₂; Item No. 14010) production in RAW 264.7 macrophages when used at concentrations ranging from 30 to 300 μ M.⁴ Avicularin (100 μ g/ml) inhibits the proliferation, migration, and invasion of, as well as induces apoptosis and cell cycle arrest at the G₀/G₁ phase in, Huh7 hepatocellular carcinoma cells.⁵

References

1. Abramson, H.N. The lipogenesis pathway as a cancer target. *J. Med. Chem.* **54**(16), 5615-5638 (2011).
2. Naeem, S., Hylands, P., and Barlow, D. Construction of an Indonesian herbal constituents database and its use in Random Forest modelling in a search for inhibitors of aldose reductase. *Bioorg. Med. Chem.* **20**(3), 1251-1258 (2012).
3. Zhang, X., Thuong, P.T., Jin, W., et al. Antioxidant activity of anthraquinones and flavonoids from flower of *Reynoutria sachalinensis*. *Arch. Pharm. Res.* **28**(1), 22-27 (2005).
4. Vo, V.A., Lee, J.-W., Chang, J.-E., et al. Avicularin inhibits lipopolysaccharide-induced inflammatory response by suppressing ERK phosphorylation in RAW 264.7 macrophages. *Biomol. Ther. (Seoul)* **20**(6), 532-537 (2012).
5. Wang, Z., Li, F., Quan, Y., et al. Avicularin ameliorates human hepatocellular carcinoma via the regulation of NF- κ B/COX-2/PPAR- γ activities. *Mol. Med. Rep.* **19**(6), 5417-5423 (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, 11/03/2022

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