

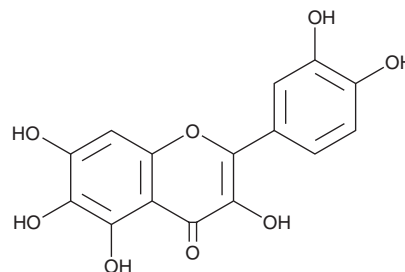
PRODUCT INFORMATION



Quercetagenin

Item No. 18645

CAS Registry No.: 90-18-6
Formal Name: 2-(3,4-dihydroxyphenyl)-3,5,6,7-tetrahydroxy-4H-1-benzopyran-4-one
Synonyms: 6-hydroxy Quercetin, NSC 115916
MF: C₁₅H₁₀O₈
FW: 318.2
Purity: ≥98%
UV/Vis.: λ_{max}: 259, 363 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

Quercetagenin is supplied as a crystalline solid. A stock solution may be made by dissolving the quercetagenin in the solvent of choice, which should be purged with an inert gas. Quercetagenin is soluble in the organic solvent DMSO.

Quercetagenin is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

The proto-oncogene serine/threonine-protein kinases, Pim-1 and Pim-2, are enzymes involved in cytokine signaling and participate in various signal transduction pathways, including cell growth, differentiation, and apoptosis. Their overexpression has been implicated in prostate cancer, some forms of leukemia, and lymphoma. Quercetagenin is a flavonol that inhibits Pim-1 with an IC₅₀ value of 0.34 μM.¹ It is selective for Pim-1, inhibiting Pim-2, PKA, RSK2, and JNK with IC₅₀ values of 3.45, 21.2, 2.82, and 4.6 μM, respectively.^{1,2} Quercetagenin has been shown to inhibit Pim-1 activity in intact RWPE2 prostate cancer cells with an ED₅₀ value of 5.5 μM, which led to significant growth inhibition.¹ It can also inhibit the growth of additional prostate epithelial cell lines at a potency proportionate to their respective level of Pim-1 protein.¹

References

1. Holder, S., Zemsanova, M., Zhang, C., *et al.* Characterization of a potent and selective small-molecule inhibitor of the PIM1 kinase. *Mol. Cancer Ther.* **6(1)**, 163-172 (2007).
2. Baek, S., Kang, N.J., Popowicz, G.M., *et al.* Structural and functional analysis of the natural JNK1 inhibitor quercetagenin. *J. Mol. Biol.* **425(2)**, 411-423 (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, 12/16/2022

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