

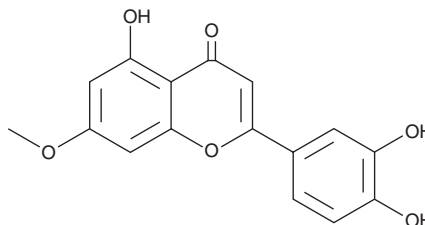
PRODUCT INFORMATION



3'-Hydroxygenkwanin

Item No. 36150

CAS Registry No.: 20243-59-8
Formal Name: 2-(3,4-dihydroxyphenyl)-5-hydroxy-7-methoxy-4H-1-benzopyran-4-one
Synonyms: HGK, Luteolin 7-methyl ether, 7-O-Methyluteolin
MF: C₁₆H₁₂O₆
FW: 300.3
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 352 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Daphne genkwa*



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

Laboratory Procedures

3'-Hydroxygenkwanin is supplied as a solid. A stock solution may be made by dissolving the 3'-hydroxygenkwanin in the solvent of choice, which should be purged with an inert gas. 3'-Hydroxygenkwanin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 3'-hydroxygenkwanin in these solvents is approximately 2 and 1 mg/ml, respectively.

Description

3'-Hydroxygenkwanin is a flavonoid that has been found in *B. balsamifera* and has antioxidant and anticancer activities.^{1,2} It scavenges superoxide anions in a cell-free assay when used at a concentration of 100 μM.¹ 3'-Hydroxygenkwanin (25 and 50 μM) decreases the levels of EGFR and phosphorylated EGFR in, and reduces viability of, PC-9 and H1975 non-small cell lung cancer (NSCLC) cells.² It also decreases intratumoral levels of EGFR and phosphorylated EGFR and reduces tumor growth in an H1975 mouse xenograft model when administered at a dose of 1 mg/kg.

References

1. Fazilatun, N., Nornisah, M., and Zhari, I. Superoxide radical scavenging properties of extracts and flavonoids isolated from the leaves of *Blumea balsamifera*. *Pharm. Biol.* **42**(6), 404–408 (2004).
2. Leu, Y.-L., Wang, T.-H., Wu, C.-C., et al. Hydroxygenkwanin suppresses non-small cell lung cancer progression by enhancing EGFR degradation. *Molecules* **25**(4), 941 (2020).

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, 10/21/2022

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