

PRODUCT INFORMATION



Recoflavone

Item No. 44582

CAS Registry No.: 203191-10-0

Formal Name: 2-[[2-(3,4-dimethoxyphenyl)-5-methoxy-4-oxo-4H-1-benzopyran-7-yl]oxy] acetic acid

Synonyms: 7-Carboxymethyloxy-3',4',5-trimethoxyflavone, DA-6034

MF: C₂₀H₁₈O₈

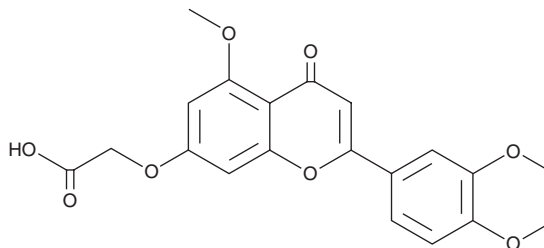
FW: 386.4

Purity: ≥98%

Supplied as: A 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

Recoflavone is supplied as a solid. A stock solution may be made by dissolving the recoflavone in the solvent of choice, which should be purged with an inert gas. Recoflavone is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution of acetonitrile:water.

Description

Recoflavone is a synthetic flavonoid with diverse biological activities and a derivative of eupatilin (Item No. 19601).¹⁻⁴ It induces chloride currents and intracellular calcium release in primary human conjunctival epithelial cells when used at concentrations of 10 and 100 μ M, respectively.¹ Recoflavone (10 and 100 μ M) induces the migration of AGS and SNU-484 human gastric epithelial cells in a scratch assay in an *in vitro* model of gastric ulcer healing.² *In vivo*, recoflavone (50 mg/kg) decreases body weight gain rate, liver weight and triglyceride levels, and hepatic macrophage and CD8⁺ T cell infiltration in a mouse model of obesity induced by a high-fat diet.³ Recoflavone (30 mg/kg) decreases disease severity, increases colon length, and increases survival in a mouse model of colitis induced by dextran sulfate (DSS; Item No. 23250).⁴ It also inhibits azoxymethane- and DSS-induced carcinogenesis in a mouse model of colitis-associated cancer.

References

1. Yang, Y.M., Park, S., Ji, H., *et al.* DA-6034 Induces [Ca²⁺]_i increase in epithelial cells. *Korean J. Physiol. Pharmacol.* **18**(2), 89-94 (2014).
2. Kim, Y.W., Lee, W.H., Choi, S.M., *et al.* DA6034 promotes gastric epithelial cell migration and wound-healing through the mTOR pathway. *J. Gastroenterol. Hepatol.* **27**(2), 397-405 (2012).
3. Kim, H.M., Kwon, M.-H., Lee, E.S., *et al.* DA-6034 ameliorates hepatic steatosis and inflammation in high fat diet-induced obese mice. *J. Yeungnam Med. Sci.* **41**(2), 103-112 (2024).
4. Nam, S.Y., Kim, J.S., Kim, J.M., *et al.* DA-6034, a derivative of flavonoid, prevents and ameliorates dextran sulfate sodium-induced colitis and inhibits colon carcinogenesis. *Exp. Biol. Med. (Maywood)* **233**(2), 180-191 (2008).

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, 11/04/2025

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