

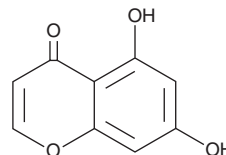
# PRODUCT INFORMATION



## 5,7-Dihydroxychromone

Item No. 29006

**CAS Registry No.:** 31721-94-5  
**Formal Name:** 5,7-dihydroxy-4H-1-benzopyran-4-one  
**MF:** C<sub>9</sub>H<sub>6</sub>O<sub>4</sub>  
**FW:** 178.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251, 258, 296 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Garcinia mangostana* L.



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

### Laboratory Procedures

5,7-Dihydroxychromone is supplied as a crystalline solid. A stock solution may be made by dissolving the 5,7-dihydroxychromone in the solvent of choice, which should be purged with an inert gas. 5,7-Dihydroxychromone is soluble in the organic solvent methanol.

### Description

5,7-Dihydroxychromone is a flavone decomposition product that has been found in peanuts and has antioxidant activities.<sup>1</sup> It inhibits increases in apoptosis and reactive oxygen species (ROS) induced by 6-OHDA (Item No. 25330) in SH-SY5Y neuroblastoma cells when used at concentrations of 0.4, 2, and 10 μM. 5,7-Dihydroxychromone (0.08-10 μM) increases the levels of HO-1, NQO1, and GCLc in SH-SY5Y cells.

### Reference

1. Kim, D.-W., Lee, K.-t., Kwon, J., *et al.* Neuroprotection against 6-OHDA-induced oxidative stress and apoptosis in SH-SY5Y cells by 5,7-Dihydroxychromone: Activation of the Nrf2/ARE pathway. *Life Sci.* **130**, 25-30 (2015).

#### 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, 08/25/2023

#### 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