

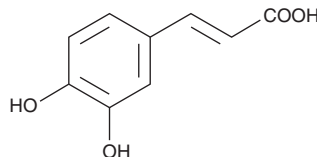
# PRODUCT INFORMATION



## Caffeic Acid

Item No. 70602

**CAS Registry No.:** 331-39-5  
**Formal Name:** 3-(3,4-dihydroxyphenyl)-2-propenoic acid  
**Synonym:** 3,4-Dihydroxycinnamic Acid  
**MF:** C<sub>9</sub>H<sub>8</sub>O<sub>4</sub>  
**FW:** 180.2  
**Purity:** ≥97%  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥2 years



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

### Laboratory Procedures

Caffeic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the caffeic acid in the solvent of choice. Caffeic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of caffeic acid in these solvents is approximately 7 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free aqueous solution is needed, it can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of caffeic acid in PBS (pH 7.2) is approximately 0.65 mg/ml. Store aqueous solutions of caffeic acid on ice and use within 12 hours of preparation. Although the aqueous solutions of caffeic acid may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

### Description

Caffeic acid inhibits both 5- and 12-lipoxygenase (LO) in a dose-dependent manner.<sup>1</sup> Depending on the type of assay, the reported IC<sub>50</sub> for 5-LO inhibition is between 3.7 μM and 72 μM.<sup>1-3</sup> The inhibition of 12-LO by caffeic acid also shows variation; its IC<sub>50</sub> has been reported between 5.1 μM and 30 μM.<sup>1-3</sup>

### References

1. Koshihara, Y., Neichi, T., Murota, S., *et al.* Caffeic acid is a selective inhibitor for leukotriene biosynthesis. *Biochim. Biophys. Acta* **792**, 92-97 (1984).
2. Kohyama, N., Nagata, T., Fujimoto, S., *et al.* Inhibition of arachidonate lipoxygenase activities by 2-(3,4-dihydroxyphenyl)ethanol, a phenolic compound from Olives. *Biosci. Biotech. Biochem.* **61**, 347-350 (1997).
3. Rao, C.V., Desai, D., Simi, B., *et al.* Inhibitory effect of caffeic acid esters on azoxymethane-induced biochemical changes and aberrant crypt foci formation in rat colon. *Cancer Res.* **53**, 4182-4188 (1993).

#### 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, 09/20/2019

#### 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