

PRODUCT INFORMATION



Chicoric Acid

Item No. 24960

CAS Registry No.: 70831-56-0
Formal Name: (2R,3R)-2,3-bis[[(2E)-3-(3,4-dihydroxyphenyl)-1-oxo-2-propen-1-yl]oxy]butanedioic acid
Synonyms: L-Chicoric Acid, Dicafeoyltartaric Acid, NSC 99173

MF: C₂₂H₁₈O₁₂

FW: 474.4

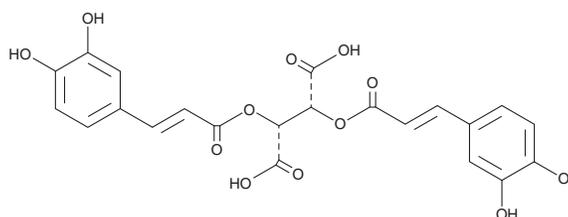
Purity: ≥95%

UV/Vis.: λ_{max}: 220, 248, 334 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

Chicoric acid is supplied as a crystalline solid. A stock solution may be made by dissolving the chicoric acid in the solvent of choice, which should be purged with an inert gas. Chicoric acid is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of chicoric acid in these solvents is approximately 25 and 30 mg/ml, respectively.

Chicoric acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, chicoric acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Chicoric acid has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Chicoric acid is a dicafeoyl ester that has been found in *C. intybus* with diverse biological activities.¹ Chicoric acid (50-200 µg/ml) dose-dependently reduces the viability of Caco-2 and HCT116 human colorectal cancer cells.² It inhibits HIV integrase activities, including 3'-processing of a DNA oligonucleotide and integration with template DNA (IC₅₀s = 1.1 and 0.8 µM, respectively).³ Chicoric acid (0.5-10 µM) noncompetitively inhibits integration of HIV DNA by HIV integrase and, at concentrations greater than or equal to 5 µM, inhibits HIV entry into H9 cells.⁴ Oral administration of chicoric acid (10 and 30 mg/kg) reduces hepatic lipid accumulation, lipid peroxidation, and fibrosis, inhibits production of pro-inflammatory cytokines and activation of NF-κB, and activates the AMPK signaling pathway in a mouse model of non-alcoholic steatohepatitis (NASH) induced by a methionine and choline-deficient diet.⁵ Chicoric acid (2 mg/kg) also reduces blood glucose levels by 54% in mice with streptozotocin-induced diabetes.⁶

References

1. Xiao, H., Xie, G., Wang, J., et al. *Food Res. Int.* **54**(1), 345-353 (2013).
2. Tsai, Y.-L., Chiu, C.-C., Chen, J.Y.-F., et al. *J. Ethnopharmacol.* **143**(3), 914-919 (2012).
3. Lin, Z., Neamati, N., Zhao, H., et al. *J. Med. Chem.* **42**(8), 1401-1414 (1999).
4. Reinke, R.A., Lee, D.J., McDougall, B.R., et al. *Virology* **326**(2), 203-219 (2004).
5. Kim, M., Yoo, G., Randy, A., et al. *Mol. Nutr. Food Res.* **61**(5), (2017).
6. Casanova, L.M., da Silva, D., Sola-Penna, M., et al. *Filoterapia* **93**, 132-141 (2014).

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, 12/06/2022

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