

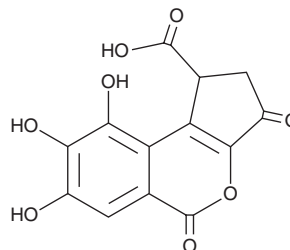
PRODUCT INFORMATION



Brevifolincarboxylic Acid

Item No. 34922

CAS Registry No.: 18490-95-4
Formal Name: 1,2,3,5-tetrahydro-7,8,9-trihydroxy-3,5-dioxo-cyclopenta[c][2]benzopyran-1-carboxylic acid
Synonym: Brevifolin Carboxylic Acid
MF: C₁₃H₈O₈
FW: 292.2
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 279, 353 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Erodii herba*



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

Laboratory Procedures

Brevifolincarboxylic acid is supplied as a solid. A stock solution may be made by dissolving the brevifolincarboxylic acid in the solvent of choice, which should be purged with an inert gas. Brevifolincarboxylic acid is soluble in DMSO.

Description

Brevifolincarboxylic acid is a phenol that has been found in *D. chrysantha* and has anticancer and anti-inflammatory activities.^{1,2} It inhibits the proliferation of PC-14 lung cancer cells (IC₅₀ = 3.95 μg/ml).¹ Brevifolincarboxylic acid (20 μM) reduces LPS-induced increases in nitric oxide (NO) production in RAW 264.7 macrophages.²

References

1. Lee, I.R. and Yang, M.Y. Phenolic compounds from *Duchesnea chrysantha* and their cytotoxic activities in human cancer cell. *Arch. Pharm. Res.* **17(6)**, 476-479 (1994).
2. Lee, B.Z., Kim, C.-S., Jeong, S.-K., et al. Anti-inflammatory isocoumarins from the bark of *Fraxinus chinensis* subsp. *rhyrachophylla*. *Nat. Prod. Res.* 1-8 (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, 12/08/2022

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