

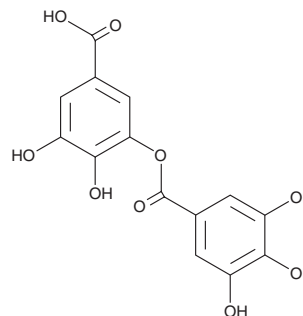
PRODUCT INFORMATION



Digallic Acid

Item No. 21224

CAS Registry No.: 536-08-3
Formal Name: 3,4-dihydroxy-5-[(3,4,5-trihydroxybenzoyl)oxy]-benzoic acid
Synonyms: *m*-Galloylgallic acid, NSC 59263
MF: C₁₄H₁₀O₉
FW: 322.2
Purity: ≥95% (mixture of *para* and *meta* isomers)
Supplied as: A solid
Storage: 4°C
Stability: ≥1 year



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

Laboratory Procedures

Digallic acid is supplied as a solid. A stock solution may be made by dissolving the digallic acid in the solvent of choice, which should be purged with an inert gas. Digallic acid is soluble in organic solvents such as methanol and DMSO.

Description

Digallic acid is a natural polyphenolic that can be produced by hydrolysis of gallotannins. Like other polyphenolic compounds, digallic acid has antioxidant activity that can be cytoprotective.¹ Digallic acid inhibits reverse transcriptases from human immunodeficiency virus ($K_i = 0.58 \mu\text{M}$) and murine leukemia virus.² It also inhibits calcium-activated chloride channels, blocking the initial agonist-stimulated chloride current.³

References

1. Bhourri, W., Derbel, S., Skandrani, I., *et al.* Study of genotoxic, antigenotoxic and antioxidant activities of the digallic acid isolated from *Pistacia lentiscus* fruits. *Toxicol. In Vitro* **24(2)**, 509-515, (2010).
2. Nakane, H., Fukushima, M., and Ono, K. Differential inhibition of reverse transcriptase and various DNA polymerases by digallic acid and its derivatives. *J. Nat. Prod.* **53(5)**, 1990 1234-1240 (1990).
3. Namkung, W., Phuan, P.W., and Verkman, A.S. TMEM16A inhibitors reveal TMEM16A as a minor component of calcium-activated chloride channel conductance in airway and intestinal epithelial cells. *J. Biol. Chem.* **286(3)**, 2365-2374 (2011).

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, 12/08/2021

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