

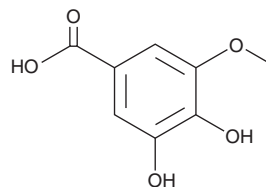
PRODUCT INFORMATION



3-O-Methylgallic Acid

Item No. 42530

CAS Registry No.: 3934-84-7
Formal Name: 3,4-dihydroxy-5-methoxy-benzoic acid
Synonyms: 4,5-Dihydroxy-3-methoxybenzoic Acid, Gallic Acid 3-methyl ester
MF: C₈H₈O₅
FW: 184.1
Purity: ≥95%
Supplied as: A 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

3-O-Methylgallic acid is supplied as a solid. A stock solution may be made by dissolving the 3-O-methylgallic acid in the solvent of choice, which should be purged with an inert gas. 3-O-Methylgallic acid is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

3-O-Methylgallic acid is a polyphenol that has been found in Keishibukuryogan-ka-yokuinin (KBGY), a traditional Japanese medicine, and has antioxidant activity.¹ It inhibits lipid hydroperoxide (LPO) generation in a cell-free assay when used at a concentration of 10 μM.

Reference

1. Matsubara, Y., Matsumoto, T., Sekiguchi, K., *et al.* Oral administration of the Japanese traditional medicine Keishibukuryogan-ka-yokuinin decreases reactive oxygen metabolites in rat plasma: Identification of chemical constituents contributing to antioxidant activity. *Molecules* **22**(2), 256 (2017).

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, 01/03/2025

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