

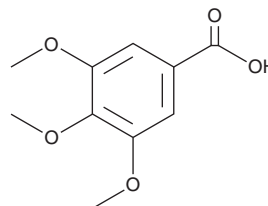
PRODUCT INFORMATION



3,4,5-Trimethoxybenzoic Acid

Item No. 34750

CAS Registry No.: 118-41-2
Formal Name: 3,4,5-trimethoxy-benzoic acid
Synonyms: Eudesmic Acid, Gallic Acid trimethyl ether, NSC 2525, Trimethylgallic Acid, TMBA
MF: C₁₀H₁₂O₅
FW: 212.2
Purity: ≥98%
UV/Vis.: λ_{max}: 214 nm
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

3,4,5-Trimethoxybenzoic acid (TMBA) is supplied as a solid. A stock solution may be made by dissolving the TMBA in the solvent of choice, which should be purged with an inert gas. TMBA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of TMBA in these solvents is approximately 30 mg/ml.

TMBA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TMBA should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. TMBA has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TMBA is an active metabolite of the opioid receptor agonist and prokinetic agent trimebutine (Item No. 31574).¹ It is active against *S. aureus* (MIC = 0.97 µg/ml).²

References

1. Wang, H., Zhou, H., Horimoto, S., *et al.* Quantitative determination of trimebutine maleate and its three metabolites in human plasma by liquid chromatography-tandem mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **779(2)**, 173-187 (2002).
2. Bisignano, G., Sanogo, R., Marino, A., *et al.* Antimicrobial activity of *Mitracarpus scaber* extract and isolated constituents. *Lett. Appl. Microbiol.* **30(2)**, 105-108 (2000).

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