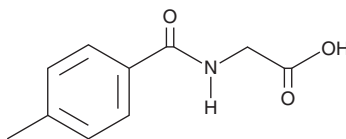


PRODUCT INFORMATION

4-Methylhippuric Acid

Item No. 46067

CAS Registry No.: 27115-50-0
Formal Name: N-(4-methylbenzoyl)-glycine
Synonyms: *para*-Methylhippuric Acid,
p-Methylhippuric Acid, 4-MHA,
 NSC 126814, *p*-Toluric Acid
MF: C₁₀H₁₁NO₃
FW: 193.2
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

4-Methylhippuric acid (4-MHA) is supplied as a solid. A stock solution may be made by dissolving the 4-MHA in the solvent of choice, which should be purged with an inert gas. 4-MHA is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 4-MHA can be prepared by directly dissolving the solid in aqueous buffers. 4-MHA is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

4-MHA is a metabolite of the volatile organic compound (VOC) and organic solvent xylene.¹ It has been used as a urinary biomarker for xylene exposure.² Higher urinary levels of 3- and 4-MHA are associated with an increased risk of high-frequency hearing loss and of osteoporosis.^{3,4}

References

1. Szűcs, S., Tóth, L., Legóza, J., *et al.* Simultaneous determination of styrene, toluene, and xylene metabolites in urine by gas chromatography/mass spectrometry. *Arch. Toxicol.* **76**(10), 560-569 (2002).
2. da Deus Honório, T.C., de Oliveira Neto, J.R., Oliveira, F.N.M., *et al.* Occupational exposure evaluation of Brazil university community to the volatile organic compounds. *J. Pharm. Biomed. Anal.* **191**, 113637 (2020).
3. Liu, Y., Long, Z., Qiu, J., *et al.* Combined effects of benzene, toluene, xylene, ethylbenzene, and styrene exposure on hearing loss mediated by oxidative stress at realistic low levels. *Environ. Pollut.* **363**(Pt 1), 125149 (2024).
4. Zhou, H., Cui, Z., Di, D., *et al.* Connecting volatile organic compounds exposure to osteoporosis risk via oxidative stress based on adverse outcome pathway methodology. *J. Environ. Sci. (China)* **155**, 806-817 (2025).

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, 07/17/2026

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