

# PRODUCT INFORMATION

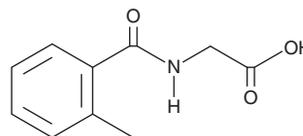


## 2-Methylhippuric Acid

Item No. 34376

**CAS Registry No.:** 42013-20-7  
**Formal Name:** N-(2-methylbenzoyl)-glycine  
**Synonyms:** 2-Methylbenzoylglycine, *o*-Methylhippuric Acid, *ortho*-Methylhippuric Acid, 2-MHA, NSC 163983, *o*-Toluric Acid

**MF:** C<sub>10</sub>H<sub>11</sub>NO<sub>3</sub>  
**FW:** 193.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 204 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

2-Methylhippuric acid is supplied as a solid. A stock solution may be made by dissolving the 2-methylhippuric acid in the solvent of choice, which should be purged with an inert gas. 2-Methylhippuric acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2-methylhippuric acid in these solvents is approximately 15, 20, and 30 mg/ml, respectively.

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 2-methylhippuric acid can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 2-methylhippuric acid in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2-Methylhippuric acid is a metabolite of the organic solvent xylene.<sup>1</sup> It has been found in the urine of industrial workers exposed to xylene.

### Reference

1. Szűcs, S., Tóth, L., Legoza, 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).

#### 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, 12/06/2022

#### 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