

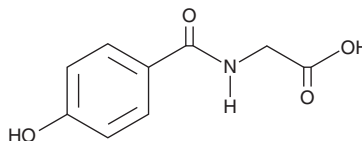
PRODUCT INFORMATION



4-Hydroxyhippuric Acid

Item No. 30115

CAS Registry No.: 2482-25-9
Formal Name: N-(4-hydroxybenzoyl)-glycine
Synonyms: *p*-Hydroxyhippuric Acid,
para-Hydroxyhippuric Acid
MF: C₉H₉NO₄
FW: 195.2
Purity: ≥98%
UV/Vis.: λ_{max}: 254 nm
Supplied as: A crystalline 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-Hydroxyhippuric acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-hydroxyhippuric acid in the solvent of choice, which should be purged with an inert gas. 4-Hydroxyhippuric acid is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 1 mg/ml.

4-Hydroxyhippuric acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-hydroxyhippuric acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 4-Hydroxyhippuric acid has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-Hydroxyhippuric acid is a dihydroxy phenolic acid inhibitor of organic anion transporter 1 (OAT1; IC₅₀ = 25 μM in opossum kidney cells).¹ It increases the production of free radicals in opossum kidney cells expressing human OAT1. 4-Hydroxyhippuric acid reduces the secretion of TNF-α, but not IL-1β or IL-6, in isolated human peripheral blood mononuclear cells (PBMCs) stimulated with LPS by 30.4% when used at a concentration of 1 μM.²

References

1. Motojima, M., Hosokawa, A., Yamato, H., *et al.* Uraemic toxins induce proximal tubular injury via organic anion transporter 1-mediated uptake. *Br. J. Pharmacol.* **135**(2), 555-563 (2002).
2. Monagas, M., Khan, N., Andrés-Lacueva, C., *et al.* Dihydroxylated phenolic acids derived from microbial metabolism reduce lipopolysaccharide-stimulated cytokine secretion by human peripheral blood mononuclear cells. *Br. J. Nutr.* **102**(2), 201-206 (2009).

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, 11/23/2022

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