

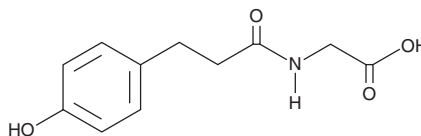
PRODUCT INFORMATION



4-Hydroxyphenylpropionylglycine

Item No. 36389

CAS Registry No.: 3850-43-9
Formal Name: N-[3-(4-hydroxyphenyl)-1-oxopropyl]-glycine
Synonyms: Phloretic Acid Glycine Conjugate, Phloretylglycine, 4-hydroxy PPG
MF: C₁₁H₁₃NO₄
FW: 223.2
Purity: ≥95%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

4-Hydroxyphenylpropionylglycine is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxyphenylpropionylglycine in the solvent of choice, which should be purged with an inert gas. 4-Hydroxyphenylpropionylglycine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-hydroxyphenylpropionylglycine in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

4-Hydroxyphenylpropionylglycine is a metabolite of the conditionally essential amino acid tyrosine.¹ It is formed from tyrosine *via* aromatic amino acid aminotransferase, tyrosine aminotransferase, and gut microbiota, followed by glycine conjugation. 4-Hydroxyphenylpropionylglycine is also a metabolite of the phenol phloretin (Item No. 14452).²

References

1. Anesi, A., Berding, K., Clarke, G., *et al.* Metabolomic workflow for the accurate and high-throughput exploration of the pathways of tryptophan, tyrosine, phenylalanine, and branched-chain amino acids in human biofluids. *J. Proteome Res.* **21**(5), 1262-1275 (2022).
2. Monge, P., Solheim, E., and Scheline, R.R. Dihydrochalcone metabolism in the rat: Phloretin. *Xenobiotica* **14**(2), 917-924 (1984).

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, 11/02/2023

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