

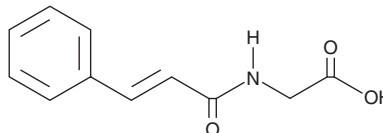
PRODUCT INFORMATION



Cinnamoylglycine

Item No. 36378

CAS Registry No.: 16534-24-0
Formal Name: N-(1-oxo-3-phenyl-2-propen-1-yl)-glycine
MF: $C_{11}H_{11}NO_3$
FW: 205.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 205, 217, 223, 274 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Cinnamoylglycine is supplied as a solid. A stock solution may be made by dissolving the cinnamoylglycine in the solvent of choice, which should be purged with an inert gas. Cinnamoylglycine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of cinnamoylglycine in these solvents is approximately 1, 25, and 15 mg/ml, respectively.

Description

Cinnamoylglycine is a metabolite of the unsaturated carboxylic acid *trans*-cinnamic acid.¹ Urinary levels of cinnamoylglycine are decreased in a mouse model of gut microbiota colonization resistance induced by piperacillin (Item No. 20766) and tazobactam (Item Nos. 25679 | 17185).² Serum levels of cinnamoylglycine are increased in patients with chronic kidney disease (CKD).³

References

1. Yan, K., Chen, W., Zhu, H., *et al.* The changes of serum metabolites in diabetic GK rats after ileal transposition surgery. *Obes. Surg.* **29**(3), 882-890 (2019).
2. Obrenovich, M.E., Tima, M., Polinkovsky, A., *et al.* Targeted metabolomics analysis identifies intestinal microbiota-derived urinary biomarkers of colonization resistance in antibiotic-treated mice. *Antimicrob. Agents Chemother.* **61**(8), e00477-e00417 (2017).
3. Boelaert, J., t'Kindt, R., Schepers, E., *et al.* State-of-the-art non-targeted metabolomics in the study of chronic kidney disease. *Metabolomics* **10**(3), 425-442 (2014).

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, 05/16/2023

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