

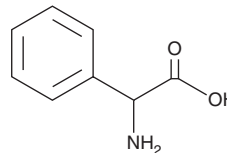
PRODUCT INFORMATION



Phenylglycine

Item No. 37186

CAS Registry No.: 2835-06-5
Formal Name: α -amino-benzeneacetic acid
Synonyms: 2-Phenylglycine, α -Phenylglycine, DL-Phenylglycine, NSC 7928, NSC 24619, NSC 32070
MF: $C_8H_9NO_2$
FW: 151.2
Purity: $\geq 95\%$
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

Phenylglycine is supplied as a solid. A stock solution may be made by dissolving the phenylglycine in the solvent of choice, which should be purged with an inert gas. Phenylglycine is soluble in 1N HCl at a concentration of approximately 10 mg/ml.

Description

Phenylglycine is a building block that has been used in the synthesis of hepatitis C virus (HCV) non-structural protein 3 (NS3) protease inhibitors and compounds with fungicidal activity.^{1,2}

References

1. Ortqvist, P., Peterson, S.D., Kerblom, E., *et al.* Phenylglycine as a novel P2 scaffold in hepatitis C virus NS3 protease inhibitors. *Bioorg. Med. Chem.* **15(3)**, 1448-1474 (2007).
2. Zhang, H., Zhao, C., Zheng, H., *et al.* Design, synthesis and bioassay of 2-phenylglycine derivatives as potential pesticide candidates. *Chem. Biodivers.* **20(1)**, e202200957 (2023).

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, 05/17/2023

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