

PRODUCT INFORMATION



H-AAA-OH

Item No. 42939

CAS Registry No.: 5874-90-8
Formal Name: L-alanyl-L-alanyl-L-alanine
Synonyms: Ala3, Ala-Ala-Ala,
L-Alanine-L-alanine-L-alanine,
Trialanine, Tri-L-alanine

Peptide Sequence: AAA-OH

MF: C₉H₁₇N₃O₄

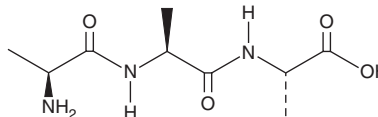
FW: 231.3

Purity: ≥95%

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

H-AAA-OH is supplied as a solid. Aqueous solutions of H-AAA-OH can be prepared by directly dissolving the solid in aqueous buffers. H-AAA-OH is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

H-AAA-OH is a tripeptide. It is a source of nitrogen for plants and microorganisms in soil.¹

Reference

1. Farrell, M.A., Hill, P.W., Farrar, J.T., *et al.* Oligopeptides represent a preferred source of organic N uptake: A global phenomenon? *Ecosystems* **16**, 133-145 (2012).

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, 04/29/2025

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