

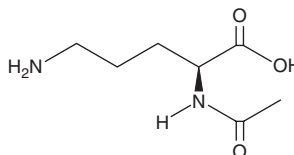
PRODUCT INFORMATION



N-acetyl-L-Ornithine

Item No. 36092

CAS Registry No.: 6205-08-9
Formal Name: N²-acetyl-L-ornithine
Synonym: N^α-acetyl-L-Ornithine
MF: C₇H₁₄N₂O₃
FW: 174.2
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

N-acetyl-L-Ornithine is supplied as a solid. A stock solution may be made by dissolving the N-acetyl-L-ornithine in the solvent of choice, which should be purged with an inert gas. N-acetyl-L-Ornithine is slightly soluble in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N-acetyl-L-ornithine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-acetyl-L-ornithine in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-acetyl-L-ornithine is an intermediate in arginine and proline metabolism and a precursor in the synthesis of the protein oxidation product γ -glutamic semialdehyde.^{1,2}

References

1. Akagawa, M., Sasaki, D., Kurota, Y., *et al.* Formation of α -amino adipic and γ -glutamic semialdehydes in proteins by the Maillard reaction. *Ann. N. Y. Acad. Sci.* **1043(1)**, 129-134 (2005).
2. Soladoye, O.P., Shand, P., Dugan, M.E.R., *et al.* Influence of cooking methods and storage time on lipid and protein oxidation and heterocyclic aromatic amines production in bacon. *Food Res. Int.* **99(1)**, 660-669 (2017).

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, 11/28/2022

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