

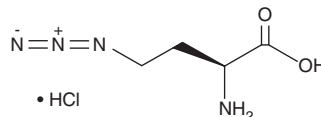
PRODUCT INFORMATION



L-Azidohomoalanine (hydrochloride)

Item No. 11786

CAS Registry No.: 942518-29-8
Formal Name: 2-amino-4-azido-butanoic acid, monohydrochloride
Synonym: L-AHA
MF: C₄H₈N₄O₂ • HCl
FW: 180.6
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

L-Azidohomoalanine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the L-azidohomoalanine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. L-Azidohomoalanine (hydrochloride) is soluble in methanol. L-Azidohomoalanine (hydrochloride) is slightly soluble in acetonitrile.

L-Azidohomoalanine (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

L-Azidohomoalanine is a clickable form of the non-proteinogenic amino acid L-homoalanine (L-2-aminobutyric acid; Item No. 38750).¹ It is composed of an L-methionine derivative functionalized with a clickable azide group. L-Azidohomoalanine has been used for semiquantitative analysis of newly synthesized proteins in A549 epithelial cancer cells and neural progenitor cells derived from patients with fragile X syndrome.^{1,2}

References

1. Wang, F., Kong, H., Meng, X., *et al.* A light-initiated chemical reporter strategy for spatiotemporal labeling of biomolecules. *RSC Chem. Biol.* **3(5)**, 539-545 (2022).
2. Raj, N., McEachin, Z.T., Harousseau, W., *et al.* Cell-type-specific profiling of human cellular models of fragile X syndrome reveal PI3K-dependent defects in translation and neurogenesis. *Cell Rep.* **35(2)**, 108991 (2021).

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, 12/18/2023

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