

4-Azido-L- homoalanine HCl (L- AHA) (100 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-AA005-100

100mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (S)-2-Amino-4-azidobutanoic acid hydrochloride

Appearance white powder

Application Notes Proteins synthesis monitoring

CAS 942518-29-8

Formula $C_4H_8N_4O_2 \cdot HCl$

MW 144.13 + 36.45

Purity > 99 % (TLC, Titration)

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com