

His(6) monoclonal antibody (6-His) (purified)

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ENZ-ABS143A-0200	200µl
------------------	-------

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 His-tag

Application IF, IP, WB

Clone 6-His

Formulation Liquid. Phosphate-buffered solution containing 0.03% thimerosal.

Host Mouse

Immunogen Recombinant protein containing the sequence HHHHHH.

Isotype IgG1

Purity Detail Affinity chromatography purified.

Recommendation Immunofluorescence (1:1,000)Immunoprecipitation (1:150)Western blot
Dilutions/Conditions (1:1,000)Optimal conditions must be determined individually for each application.

Specificity Recognizes proteins containing the HHHHHH sequence. The sequence of the protein flanking the poly His-tag does not influence antibody binding.

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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