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.

lsotype lgG1

Purity Detail Affinity chromatography purified.

Recommendation Dilutions/Conditions Immunofluorescence (1:1,000)Immunoprecipitation

(1:150)Western blot (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

•

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com