

His(6) Recombinant monoclonal antibody (Hamster IgGκ)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Polyhistidine tag
Application	ELISA, WB
Clone	3D5
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Hamster (Armenian)
Immunogen	Purified his tagged Fab fragment.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Specificity	Recognizes carboxy-terminal oligo-histidine tags (6x His tags) of fusion proteins. It also has affinity for internal His-tags but no affinity for N-terminal His-tags.
Technical Info / Product Notes	This is a chimeric hamster antibody, based on the original mouse IgG2b format, created for improved compatibility with existing reagents, assays and techniques.
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
[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)