

His(6) Recombinant monoclonal antibody (Mouse IgG2bk)

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

Ordering Information

[Order Online »](#)

ENZ-ABS535-0200	200µg
-----------------	-------

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	Mouse
Immunogen	Purified his tagged Fab fragment.
Isotype	IgG2bk
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.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee