His(6) Recombinant monoclonal antibody (Human IgG3κ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS539-0200

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 Human

Immunogen Purified his tagged Fab fragment.

lgG3κ

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 NotesThis is a chimeric human 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



info-

eu@enzolifesciences.com