

Anti-His

Mouse Monoclonal Antibody

Catalog # H99-61M

Lot # Y970-3

Cited Applications

- Western blot (1:1000)
- Immunostaining
- ELISA
- Immunoprecipitation

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the N-terminal, C-terminal or internal 6X His-tagged fusion protein

Cross Reactivity

- All proteins tagged with 6X His

Host/Isotype/Clone#

Mouse, IgG2b, EH158

Immunogen

HHHHHH peptide conjugated to KLH

Formulation

10mM PBS pH 7.2

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

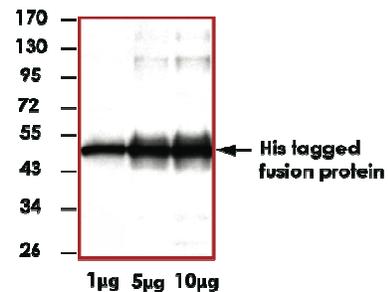
Scientific Background

This polyhistidine epitope tag is generally comprised of six consecutive histidine amino acid residues located at the N-terminal, C-terminal, or internally [1]. The 6 His-Tag is widely used because of its affinity to bind nickel or cobalt metal ions attached to sepharose, which can then be used to purify the protein in a native or denatured state [2].

References

1. Jonathan, W. et al: Epitope tagging, Annual Review of Genetics, 1998; 32, 601-618.
2. Schmeisser, H, et al; Binding Characteristics of IFN-alpha Subvariants to IFNAR2-EC and Influence of the 6-Histidine Tag. J Interferon Cytokine Res. 2006; Dec. 26(12):866-76.

Sample Data



Representative western blot with Anti-His (1:1000) using 1 µg, 5 µg, 10 µg of HEK 293 cell lysate transfected with a 52 kDa His tagged fusion protein.

Anti-His

Mouse Monoclonal Antibody

Catalog # H99-61M

Lot # Y970-3

Purification Affinity purified using Protein A

Concentration 1.0 µg/µL

Stability 1 yr at -20°C from date of shipment

Storage & Shipping Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.