

Datasheet for Mouse anti His-Tag mAb

Catalog number: mAb-02620-020

Background: Consecutive histidine residues (usually 6 to 10 in length) are often inserted into the amino acid sequences of recombinant proteins. The resulting His-tagged proteins can be detected by using His antibodies.

■ Product Information

Symbol: His
Alias: His;His tag;His-tag
Cross reactivity: Species independent
Expected Protein Weight:
Speices:
Source: Mouse
Isotype: IgG1
Purity: Affinity purification
Storage: Store at -20°C. Avoid freeze / thaw cycles.|Buffer: PBS with 0.09% Sodium azide,50% glycerol,pH7.3.

■ Immunogen Information

Gene ID:
Swiss:
Immunogen: A synthetic peptide corresponding to His tag.

■ Recommended Dilutions

WB 1:5000 - 1:20000
IF/ICC 1:50-1:200
IP 0.5ug-4ug antibody for 200ug-400ug extracts of whole cells
ELISA Recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.

Citation of product: If use of this item results in a publication, please use this information: Mouse anti His-Tag mAb(mAb-02620-020; Genecopoeia, Inc, Rockville, MD).

GeneCopoeia Products are for Research Use Only
9620 Medical Center Drive, #101 Rockville, MD 20850 USA
Tel: 301-762-0888;Fax:301-762-3888
Email: support@genecopoeia.com; Web:www.genecopoeia.com

Copyright © 2024 GeneCopoeia, Inc.