

## 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](mailto:support@genecopoeia.com); Web:[www.genecopoeia.com](http://www.genecopoeia.com)

Copyright © 2024 GeneCopoeia, Inc.