

**OmicsLink™ Anti-His Tag Antibody**  
Mouse Monoclonal IgG2b

**Catalog Number**

CGAB-HIS-0100 (100 µg)  
CGAB-HIS-0050 (50 µg)

**Description**

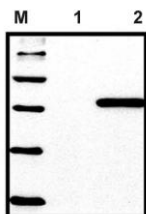
Anti-6XHis tag antibody is a purified mouse IgG2b monoclonal antibody that binds to 6XHis C- or N-terminally-tagged fusion proteins. Anti-6XHis tag antibody can be used to identify and capture fusion proteins by western blots and immunoprecipitation. Immobilized anti-6XHis is also used for affinity purification of His-tagged proteins. OmicsLink™ anti-tag antibodies are ideal for use with OmicsLink™ expression-ready ORF cDNA clones.

**Immunogen**

The OmicsLink™ anti-His tag antibody was produced by immunizing mice with a 6XHis synthetic peptide.

**Specificity**

The anti-His tag monoclonal antibody recognizes both N- or C-terminal His-tagged proteins.



**Figure 1.** 6XHis western blot analysis using 293 cells transfected with His-tagged vector (lane 2) and untransfected control (lane 1).

**Formulation**

0.25 mg/ml in PBS, pH 7.4 with 0.05% sodium azide and 50% (v/v) glycerol.

**Applications**

Western blot, immunoprecipitation, immunostaining and affinity purification.

**Storage/Stability**

Stable for 1 year at -20°C from shipment date. Briefly centrifuge the original vial prior to opening the cap to collect the product at the bottom of the tube.

**Starting dilution recommendations**

Optimal working conditions should be determined by the user. General recommendations are western blot at 1:1000; immunoprecipitation and immunostaining at 1:200

GeneCopoeia, Inc.  
9620 Medical Center Drive, #101  
Rockville, Maryland 20850  
Tel: 301-762-0888 Fax: 301-762-8333  
Email: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)  
Web: [www.genecopoeia.com](http://www.genecopoeia.com)