

OmicsLink™ Anti-HA Tag Antibody
Mouse Monoclonal IgG3

Catalog Number

CGAB-HA-0100 (100 µg)

CGAB-HA-0050 (50 µg)

Description

Anti-HA tag antibody is a purified mouse IgG3 monoclonal antibody that binds to HA-fusion proteins. Anti-HA tag antibody can be used to identify and capture fusion proteins by western blots and immunoprecipitation. Immobilized anti-HA is also used for affinity purification of HA fusion proteins. OmicsLink™ anti-tag antibodies are ideal for use with OmicsLink™ expression-ready ORF cDNA clones.

Immunogen

The OmicsLink™ anti-HA tag antibody was produced by immunizing mice with a synthetic peptide from influenza hemagglutinin epitope (YPYDVPDYA) coupled to the KLH.

Specificity

The anti-HA tag monoclonal antibody recognizes HA-tagged proteins over-expressed in cells including both N- or C-termini of targeted proteins in transfected mammalian cells.

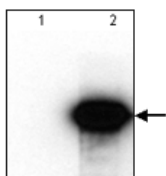


Figure 1. HA western blot analysis of 293 cells transfected with HA-tagged vector (lane 2). The untransfected control is in lane 1.

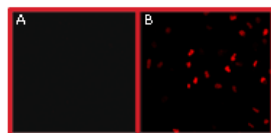


Figure 2. Immunofluorescence staining of HA-tag fusion protein in a stable expressing cell line (B) and a control cell line (A).

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
Web: www.geneCopoeia.com