

OmicsLink™ Anti-Myc Tag Antibody
Mouse Monoclonal IgG1

Catalog Number

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

Description

Anti-Myc tag is a purified mouse IgG1 monoclonal antibody that binds to cMyc-fusion proteins. Anti-Myc tag antibody can be used to identify and capture fusion proteins by western blots and immunoprecipitation. Immobilized anti-Myc is also used for affinity purification of cMyc fusion proteins. OmicsLink™ anti-tag antibodies are ideal for use with OmicsLink™ expression-ready ORF cDNA clones.

Immunogen

The OmicsLink™ anti-Myc tag antibody was produced by immunizing mice with a synthetic peptide (EQKLISEEDL) coupled to KLH.

Specificity

The anti-Myc tag monoclonal antibody recognizes over-expressed proteins containing Myc epitope tag fused to either N- or C-terminal of targeted proteins in transfected mammalian cells.

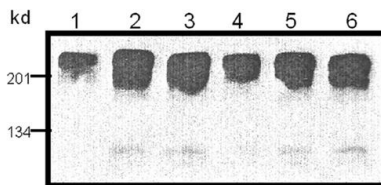


Figure 1. Western blot analysis of Myc-tagged fusion construct using different transfection reagents and variable amounts of fusion DNA construct.

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