

OmicsLink™ Anti-GFP Tag Antibody
Mouse Monoclonal IgG1

Catalog Number

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

Description

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

Immunogen

The OmicsLink™ anti-GFP tag antibody was produced by immunizing mice with a synthetic peptide of Green Fluorescent Protein (GFP) from the jellyfish *Aequorea victoria*.

Specificity

The anti-GFP tag monoclonal antibody recognizes GFP and its variants including eGFP, YFP, eYFP, and CFP.

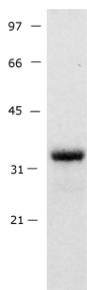


Figure 1. GFP western blot analysis of 293 cells transfected with a GFP construct. The antibody shows a 33 kDa band.

Formulation

0.25 mg/ml in PBS, pH 7.2 with 0.1% 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