

OmicsLink™ Anti-GST Tag Antibody
Mouse Monoclonal IgG2a

Catalog Number

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

Description

Anti-GST (glutathione-S-transferase) tag antibody is a purified mouse IgG2a monoclonal antibody that binds to GST-fusion proteins. Anti-GST tag antibody can be used to identify and capture fusion proteins by western blots and immunoprecipitation. Immobilized anti-GST is also used for affinity purification of GST fusion proteins. OmicsLink™ anti-tag antibodies are ideal for use with OmicsLink™ expression-ready ORF cDNA clones.

Immunogen

The OmicsLink™ anti-GST tag antibody was produced by immunizing mice with a synthetic peptide of *Schistosoma japonicum* GST.

Specificity

The anti-GST tag monoclonal antibody detects over-expressed GST fusion proteins.

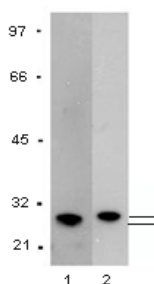


Figure 1. 20 ng of recombinant GST protein (lane 1) and 20 µg of 293 cell lysates were transfected with a GST expression vector and probed with OmicsLink Anti-GST (1 µg/ml).

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

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