

OmicsLink™ Anti-D Tag Antibody
Mouse Monoclonal IgG2b

Catalog number

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

Description

Anti-D tag antibody is a purified mouse IgG2b monoclonal antibody that binds to D epitope (DYKDDDDK, also known as FLAG®)-tagged fusion proteins. Anti-D tag antibody can be used to identify and capture fusion proteins by western blots and immunoprecipitation. Immobilized anti-D is also used for affinity purification of D-tag fusion proteins. OmicsLink™ anti-tag antibodies are ideal for use with OmicsLink™ expression-ready ORF cDNA clones.

Immunogen

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

Specificity

The anti-D tag antibody recognizes over-expressed recombinant proteins containing the D epitope tag (DYKDDDDK) fused to either the N- or C-terminal of targeted proteins in transfected mammalian cells or other expression systems.



Figure 1. Western blot analysis using 293 cells transfected with D tag vectors. Lane1: highest concentration at 0.5 mg/ml Anti-D tag antibody, Lane 4: lowest concentration at 0.1 µg/ml.

Formulation

0.25 mg/ml in PBS, pH 7.2 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 after 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