

OmicsLink™ Anti-Tag Antibodies

Overview

GeneCopoeia OmicsLink™ Anti-Tag Antibodies are monoclonal mouse IgG antibodies that bind to 6xHis-, GFP-, mCherry-, GST-, D*-, HA-, or Myc-tagged fusion proteins. The antibodies can be used to identify and capture fusion proteins by western blots, immunostaining and immunoprecipitation. The immobilized antibodies can also be used for affinity purification of tagged fusion proteins. They are ideal for use with OmicsLink™ expression-ready ORF cDNA clones, which have a wide-range of fusion tag collections.

The OmicsLink anti-tag antibodies are produced by immunizing mice with specific synthetic peptides of the fusion tags. They recognize over-expressed recombinant proteins with the epitope tags fused to either the N- or C-terminal of the protein in transfected mammalian cells or other expression systems.

Catalog number	Product	Type
CGAB-DDK-0050 CGAB-DDK-0100	OmicsLink™ Anti-D* Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG
CGAB-GFP-0050 CGAB-GFP-0100	OmicsLink™ Anti-GFP Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG
CGAB-RFP-0050 CGAB-RFP-0100	OmicsLink™ Anti-mCherry Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG
CGAB-GST-0050 CGAB-GST-0100	OmicsLink™ Anti-GST Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG
CGAB-MYC-0050 CGAB-MYC-0100	OmicsLink™ Anti-MYC Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG
CGAB-HA-0050 CGAB-HA-0100	OmicsLink™ Anti-HA Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG
CGAB-HIS-0050 CGAB-HIS-0100	OmicsLink™ Anti-HIS Antibody (50 µg or 100 µg)	Monoclonal Mouse IgG

*D tag is also known as FLAG® tag.

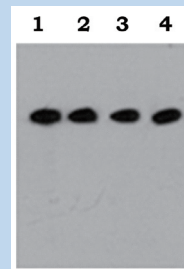


Figure 1. Western blot analysis using 293 cells transfected with D tag vectors. Lane 1: highest concentration at 0.5 mg/ml Anti-D tag antibody. Lane 4: lowest concentration at 0.1 µg/ml.

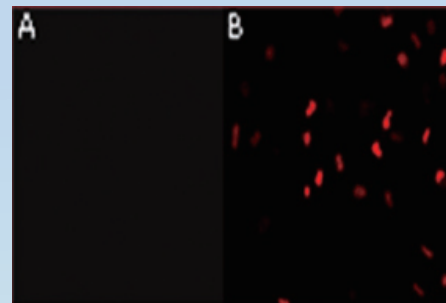
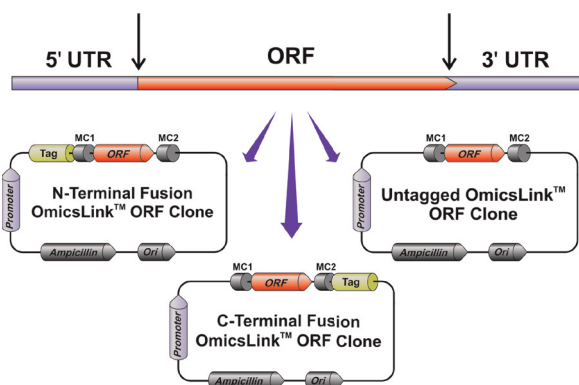


Figure 2. Immunofluorescence staining of HA-tag fusion protein in a stably expressing cell line (B) and a control cell line (A).

OmicsLink™ Anti-Tag Antibodies

Related products

OmicsLink™ expression-ready ORF cDNA clones



Expression system	Mammalian, bacterial, insect, lentiviral, yeast, wheat germ cell free
Promoter	CMV, T7, Tac, EF1 α , GAL1, pADH, AcMNPV polyhedrin, custom promoter
Selection marker	Neomycin, puromycin, hygromycin, blasticidin, zeocin
Fusion tag	<ul style="list-style-type: none"> Fluorescent tags: eGFP, eYFP, eCFP, mCherry Multifunctional tags: HaloTag®, AviTag™ Solubility and purification tags: 6xHis, SUMO, Flag, GST, MBP, 3xFlag Antibody immunoprecipitation tags: 3xHA, Myc, Flag, 3xFlag IRES- coexpressed tags: Avi+IRES-Biotin ligase, Myc+IRES-eGFP, IRES-eGFP, IRES-Neomycin, IRES-Luciferase, etc.
Vector type	Lentiviral and non-viral vectors

To order

Please visit www.genecopoeia.com or contact us directly.

GeneCopoeia, Inc.
 9620 Medical Center Drive, Suite 101
 Rockville, MD 20850, USA
 Email inquiry@genecopoeia.com
 Tel +1 (301) 762-0888
 Toll free +1 (866) 360-9531
 Fax +1 (301) 762-3888
 Website www.genecopoeia.com