

## OmicsLink™ Anti-mCherry Antibody

Mouse Monoclonal IgG

#### **Catalog Number**

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

#### **Description**

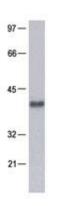
Anti-mCherry antibody is a purified mouse IgG monoclonal antibody that binds to mCherry protein. Anti-mCherry antibody can be used to identify and capture mCherry or fusion proteins by western blots and immunoprecipitation. Immobilized anti-mCherry can also be used for affinity purification of mCherry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

#### **Immunogen**

The OmicsLink<sup>™</sup> anti-mCherry antibody was raised against a purified Red Fluorescent Protein (RFP, also known as DsRed) of *Discosoma sp.* origin.

## Specificity

The anti-mCherry monoclonal antibody recognizes RFP (DsRed) and its variants including mCherry, plum, and probably other RFP-derived fluorescent proteins.



**Figure 1.** mCherry western blot analysis of 293 cells transfected with a fusion construct which contains a mCherry tag.

## **Formulation**

# 0.25 mg/ml in PBS, pH 7.2 with 0.1% sodium azide and 50% (v/v) glycerol.

# **Applications**

Western blot, ELISA, 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 1:1,000; immunoprecipitation at 5 µg/mg lysate and immunostaining at 1:200.

GeneCopoeia, Inc. 9620 Medical Center Drive, #101 Rockville, Maryland 20850

Tel: 301-762-0888 Fax: 301-762-8333 Email: <a href="mailto:inquiry@genecopoeia.com">inquiry@genecopoeia.com</a> Web: <a href="mailto:www.genecopoeia.com">www.genecopoeia.com</a>