

---

## Goat Anti-Mouse Antibody, Biotin Conjugated

(Catalog No. A12000)

---

### Description

Goat anti-mouse IgG recognizes mouse IgG whole molecule. This secondary antibody was purified using antigen affinity chromatography. The antibody is conjugated with biotin.

### Antibody Type

Polyclonal Antibody

### Purification

Liquid; this product was prepared from monospecific antiserum by immunoaffinity chromatography, followed by solid phase adsorption(s) to remove any unwanted reactivities.

### Immunogen

Mouse IgG whole molecule

### Isotype

IgG

### Size

350 µg

### Formulation

In 10 mM sodium phosphate, 75 mM NaCl, 50% (v/v) glycerol, pH 7.2 (0.01% (w/v) sodium azide, 10 mg/mL BSA

### Specificity

Mouse

### Storage

Store at 4°C. For long term storage store at -20°C. Aliquot to avoid repeated freezing and thawing. For maximum recovery of the product, centrifuge the vial prior to opening the cap.

### Handling Recommendations

The optimal working dilution should be determined by the end user.

### Applications & Suggested Dilutions

Western Blot: 1:1000-1: 10000; Immunohistochemistry: 1:100-1:500; Immunofluorescence: 1:100-1:500; ELISA: 1:2000-1:20000.

### Ordering Information:

Products	Size	Cat. No.
Goat Anti-Mouse Antibody, Biotin Conjugated	350 µg	A12000-1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*