

## Biotinylated goat anti-mouse IgG (H+L)

Catalog Number: 16729

Unit Size: 1 mg

### Product Details

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 2 mg/mL BSA

### Unit Details

---

Reconstitution Volume	1 mL ddH <sub>2</sub> O
-----------------------	-------------------------

### Antibody Properties

---

Species Reactivity	Mouse
Class	Secondary
Clonality	Polyclonal
Host	Goat

### Chemical Properties

---

Molecular Weight	~150000
------------------	---------

### Biological Properties

---

Stabilizer	2 mg/mL BSA
Appearance	Solid
Preparation	Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG. The antibody is conjugated with Biotinylated under optimal conditions.
Application	Flow Cytometry (FACS), IF, IHC, ELISA, WB
Recommended Dilutions	Suggested dilutions are only guidelines; users should titrate the product for their specific assay using appropriate controls

#### Application

#### Recommended dilution

Flow Cytometry (FACS)

1-5 µg/mL

IF	2 µg/mL
IHC	1-10 µg/mL
ELISA	100 ng/mL
WB	1-10 µg/mL

## Spectral Properties

---

Conjugate                      Biotinylated

## Applications

---

This biotinylated goat anti-mouse IgG (H+L) is affinity-purified to have high affinity. It recognizes both heavy and light chains (H+L). It is biotinylated to ensure the maximum degree of labeling without compromising the specificity or affinity of the antibody. It can be used for tissue and cell staining, in situ hybridization, Elispot ELISAs and blots.