

## MegaWox™ polyHRP-Goat Anti-Rabbit IgG Conjugate

Catalog Number: 11037

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), 0.01% thimerosal, 0.2% (w/v) BSA

### Antibody Properties

---

Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat
Immunogen	IgG

### Biological Properties

---

Appearance	Liquid
Preparation	Goat anti-Rabbit IgG (H+L) is produced in goat with pooled total Rabbit IgG, and affinity purified with Rabbit IgG coupled beads. The antibody is conjugated with MegaWox™ polyHRP under optimal condition.
Soluble In	Water
Application	IF, FACS

### Spectral Properties

---

Conjugate	polyHRP
-----------	---------

### Applications

---

Goat anti-rabbit secondary antibodies are affinity-purified antibodies with well-characterized specificity for rabbit immunoglobulins and are useful in the detection, sorting or purification of its specified target. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies can bind to a single primary antibody. MegaWox™ polyHRP-Goat Anti-Rabbit IgG Conjugate is designed to deliver the highest sensitivity and low background in immunoassays where sample volume is limited or when the target molecule is present at low levels. The goat Anti-Rabbit IgG poly-HRP

conjugate is purified to remove unconjugated goat Anti-Rabbit IgG molecules that competes for binding sites with its HRP-conjugates. In addition, the conjugate is devoid of unconjugated HRP that can cause background signal. MegaWox™ polyHRP-Goat Anti-Rabbit IgG is compatible with chromogenic, fluorogenic and chemiluminescent HRP substrates used in ELISA, Western blotting, immunohistochemistry (IHC) and nucleic acid hybridization assays. It has been validated to be used with our TSA and Styramide™ fluorescent HRP substrates for ultrasensitive detection of low abundant biological targets.