

## HRP Goat anti-Mouse IgG

**Catalog No:** 15014

**RRID:** AB\_3216396

**Application(s):** ELISA, IHC, WB

**Quantity:** 400 µg

**Purification:** Affinity Purified

**Host:** Goat

**Background:** This high quality anti-mouse IgG secondary antibody was raised in goat and has been affinity purified and conjugated to horseradish peroxidase (HRP). By improving this procedure we are able to offer antibodies with high specificity which will reduce the background and increase the sensitivity of detection in your assay.

**Immunogen:** Mouse IgG

**Buffer:** 0.01 M sodium phosphate, 0.25 M sodium chloride pH 7.6, 30% Glycerol and 15 mg/ml IgG-Free stabilizer.

**Application Notes:**

**Working dilutions:**

IHC: 1:500-1:5,000

ELISA & WB: 1:5,000-1:50,000

**Storage and Guarantee:** Avoid repeated freeze/thaw cycles by aliquoting into single-use fractions for storage at -70°C or below. Prepare working dilution on day of use. Product is stable for about 6 weeks at 2°-8°C as an undiluted liquid. Do not use sodium azide with HRP-conjugated antibodies as it will inactivate the horseradish peroxidase. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

