

# Goat Anti-Human Kappa - HRP

$\kappa$  chain specific- Revelation antibody

Catalog # hrp-igak

For research use only, not for diagnostic or therapeutic use

Version # 08D25-MT

## PRODUCT INFORMATION

### Content

1 ml of goat anti-human kappa - HRP (horseradish peroxidase conjugate) in 50% glycerol/50% PBS, pH 7.4. No preservative added.

### Storage

Product is shipped at room temperature. Store at 2-8°C or long-term at -20°C.

## DESCRIPTION

### Source

Isolated from antisera of goats hyperimmunized with human myeloma proteins containing  $\kappa$  light chains

### Cross Absorption

Pooled human myeloma proteins with  $\lambda$  light chains

### Purification

Affinity chromatography on pooled human Igs with  $\kappa$  light chains covalently linked to agarose

### Specificity

Reacts with human  $\kappa$  light chains as demonstrated by ELISA

## CHARACTERIZATION

To insure lot-to-lot consistency, each batch of product is tested by ELISA conform to characteristics of a standard reference reagent.

## APPLICATIONS

- Enzyme-Linked-Immunesorbent-Assay (ELISA)
- Western blotting
- Dot- and slot-immunoblotting

## WORKING DILUTIONS

**ELISA:** 1:4,000 - 1:8,000

**Other applications:** Since applications vary, you should determine the optimum working dilution of the product that is appropriate for your specific need.

---

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873  
Outside US: (+1) 858-457-5873  
Europe: +33 562-71-69-39  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)  
Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite 100  
San Diego, CA 92121 - USA