

Optimize your assay with solutions from ICT.

# HRP AffiPure Secondary Antibodies

## • HRP Goat anti-Rabbit IgG Fc, Catalog #6293

### Vial Composition

Each 1.0 mL vial contains 1.0 mg Goat anti-Rabbit IgG Fc conjugated to 2-4 HRP enzymes per IgG molecule. The conjugate is dissolved in a Tris-based conjugate stabilization buffer containing 50 ppm ProClin® 300 anti-microbial agent.

### Antibody Production Summary

Goats were immunized with a purified preparation of rabbit IgG Fc fragment that was obtained from a purified normal rabbit IgG antibody protein pool.

The hyperimmune IgG fraction was initially solid-phase-adsorbed to obtain class specificity. Subsequently, the anti-Fc antibody was adsorbed to and eluted from a second solid-phase immunoaffinity purification gel containing covalently coupled rabbit IgG Fc protein.

### Horseradish Peroxidase (HRP)

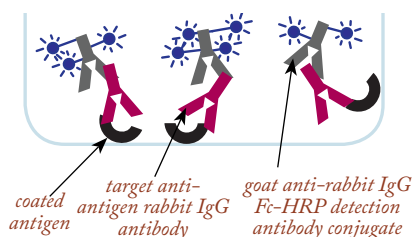
#### Conjugation Summary

Affinity purified, anti-rabbit IgG Fc specific goat IgG was covalently coupled to HRP using a maleimide (M)-facilitated conjugation technique. In this process, free sulfhydryl groups are added to the IgG Fc prep just prior to its reaction with a 4-fold molar excess of HRP-M. This process produces an IgG-HRP conjugate that contains between 2 and 4 HRP molecules per IgG molecule.

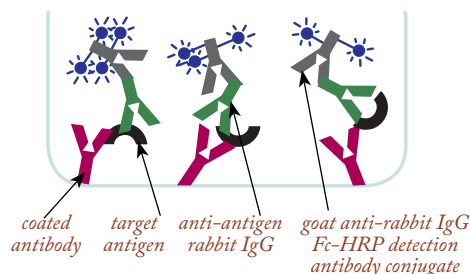
### Specificity Summary

Using ELISA and IEP analysis techniques, the IgG component of this conjugate reacts only with IgG in normal rabbit serum and does not react with other rabbit serum proteins.

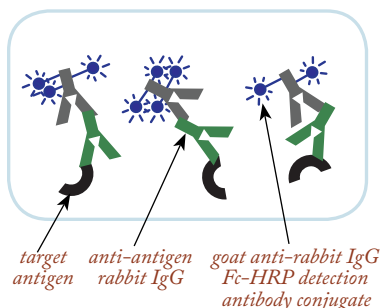
#### Antigen-Down ELISA



#### Antibody-Sandwich ELISA



#### Western Blot



### HRP Goat anti-Rabbit IgG Fc

Volume	Size	Catalog#
1 mL	1 mg	6293

### RECOMMENDED DILUTIONS

Optimal conjugate dilutions should be determined for your particular assay conditions. Here are some guidelines:

- **Western Blotting**  
1:500 – 1:2,000
- **ELISA**  
1:5,000 – 1:20,000

### SPECIFICATIONS

- Molecular composition: IgG-(HRP)2-4
- Molecular Weight: 230Kd – 310Kd
- Clear Tris buffer with 50 ppm ProClin® 300.
- Ready to dilute.

### STORAGE

- 24 months from receipt when stored at 2-8°C.

### SAFETY

- See Safety Data Sheet (SDS) for any warnings.
- SDS available at [www.immunochemistry.com](http://www.immunochemistry.com) or by calling ICT at 952-888-8788 or 1-800-829-3194.
- For research use only.
- Not for use in diagnostic procedures.

Check out more ELISA solutions on our website [www.immunochemistry.com](http://www.immunochemistry.com)

