

Optimize your assay with solutions from ICT.

# HRP AffiPure Secondary Antibodies

## • HRP Goat anti-Human IgG Fc, Catalog #6291

### Vial Composition

Each 1.0 mL vial contains 1.0 mg Goat anti-Human 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 human IgG Fc fragment that was obtained from a purified normal human 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 human IgG Fc protein.

### Horseradish Peroxidase (HRP)

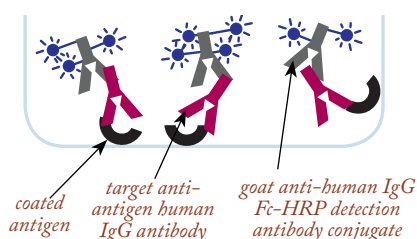
#### Conjugation Summary

Affinity purified, anti-human 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.

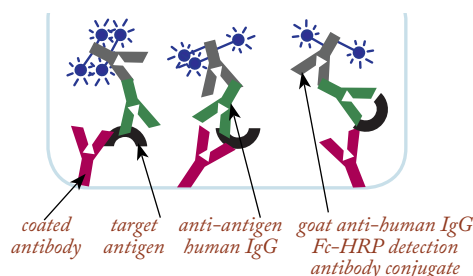
#### Specificity Summary

Using ID, IEP, and ELISA analysis techniques, the IgG component of this conjugate reacts only with the Fc fragment of IgG. It does not react with kappa or lambda light chains.

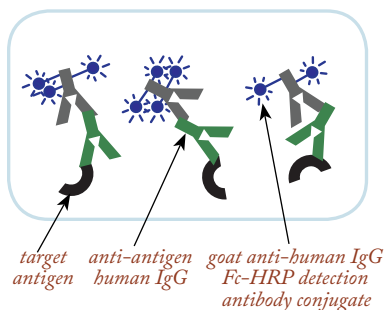
#### Antigen-Down ELISA



#### Antibody-Sandwich ELISA



#### Western Blot



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



### HRP Goat anti-Human IgG Fc

Volume	Size	Catalog#
1 mL	1 mg	6291

### 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.
- Non-reactive with k or l chains.
- 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.

For more detailed protocols, please visit  
[www.immunochemistry.com](http://www.immunochemistry.com).