

**XFD350 goat anti-rabbit IgG (H+L) \*Cross Adsorbed\***

Catalog Number: 16395

Unit Size: 1 mg

**Product Details**

---

|                    |                                            |
|--------------------|--------------------------------------------|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date    | 12 months upon receiving                   |

**Chemical Properties**

---

|                  |         |
|------------------|---------|
| Appearance       | Solid   |
| Molecular Weight | ~150000 |
| Soluble In       | Water   |

**Spectral Properties**

---

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 343 nm |
| Emission Wavelength   | 441 nm |

**Applications**

---

XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). Our goat anti-rabbit IgG whole antibodies have been cross-adsorbed against human IgG and human serum prior to conjugation to minimize cross-reactivity. This XFD350 labeled-goat anti-rabbit IgG conjugate is prepared by the reaction of cross-adsorbed goat anti-rabbit IgG whole antibody with XFD350 NHS ester, which has the same molecule to Alexa Fluor® 350 NHS ester. Each conjugate has typically 4-6 fluorophores per IgG molecule. Fluorescent secondary antibody conjugates are useful in the detection, sorting, or purification of its specified target and ideal for fluorescence microscopy and confocal laser scanning microscopy, flow cytometry, and fluorescent western detection.