

## Rabbit Anti-Cy5 Antibody, iFluor® 350-Labeled

Catalog Number: 46912

Unit Size: 100 ug

### Product Details

---

|                    |                                                |
|--------------------|------------------------------------------------|
| Storage Conditions | Refrigerated (2-8 °C), Minimize light exposure |
| Expiration Date    | 12 months upon receiving                       |
| Concentration      | 1 mg/mL                                        |
| Formulation        | PBS, BSA                                       |

### Unit Details

---

|                       |              |
|-----------------------|--------------|
| Reconstitution Volume | 0.1 mL ddH2O |
|-----------------------|--------------|

### Antibody Properties

---

|           |                                |
|-----------|--------------------------------|
| Clonality | Polyclonal                     |
| Host      | Rabbit                         |
| Immunogen | Cy5 coupled to carrier protein |

### Biological Properties

---

|             |                                                                                                                                                                             |
|-------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Stabilizer  | BSA                                                                                                                                                                         |
| Appearance  | Solid                                                                                                                                                                       |
| Preparation | It is produced with Cy5 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 350 under optimal conditions. |
| Soluble In  | Water                                                                                                                                                                       |
| Application | ELISA, IF, FACS                                                                                                                                                             |

### Spectral Properties

---

|                       |            |
|-----------------------|------------|
| Conjugate             | iFluor 350 |
| Excitation Wavelength | 345 nm     |
| Emission Wavelength   | 450 nm     |

### Applications

---

Rabbit Anti-Cy5 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Cy5-conjugated molecules and is directly labeled with iFluor® 350, a UV-excitable fluorophore with excitation/emission maxima of ~345/450 nm. This direct labeling format supports one-step detection of Cy5-labeled biomolecules without the need for a secondary antibody. The antibody may exhibit cross-reactivity with related cyanine-based dyes. It is recommended for dye conjugation validation, signal amplification in low-abundance targets, and multicolor fluorescence applications in microscopy, and flow cytometry.