

## Rabbit Anti-Cy3 Antibody, iFluor® 555-Labeled

Catalog Number: 46891

Unit Size: 100 ug

### Product Details

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.2% (w/v) BSA

### Unit Details

---

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

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Cy3 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	Antibody is produced with immunogen in rabbit, affinity purified and conjugated with iFluor® 555 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

### Applications

---

Rabbit Anti-Cy3 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Cy3-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore excited at around 550 nm. The direct labeling format enables one-step fluorescence detection, reducing background and improving workflow efficiency.

The antibody may cross-react with related dyes such as Alexa Fluor® 555, ATTO 550, and iFluor® 555. Its spectral alignment with Cy3 allows enhanced signal amplification without additional detection reagents, making it suitable for multiplex assays, microscopy, and flow cytometry.

*Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.*