

## Rabbit Anti-Cy3 Antibody, Biotinylated

Catalog Number: 46900

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 ddH <sub>2</sub> O
-----------------------	---------------------------

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Cy3 coupled to carrier protein

### Chemical Properties

---

Molecular Weight	~150000
------------------	---------

### Biological Properties

---

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

### Applications

---

Rabbit Anti-Cy3 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Cy3-labeled targets and is directly conjugated to biotin. This biotinylated format enables flexible signal detection using streptavidin- or avidin-linked reporters across a wide range of fluorescence and enzymatic assays.

The antibody may cross-react with structurally related fluorophores such as Alexa Fluor® 555, iFluor® 555, and ATTO 550. It is ideal for signal amplification in low-abundance target detection, verifying Cy3 conjugation efficiency, and performing multiplex fluorescence workflows.

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