

## Rabbit Anti-Alexa Fluor 555 Antibody, Biotinylated

Catalog Number: 46550

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	0.1 mL ddH <sub>2</sub> O
Volume	

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 555 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 555 coupled carrier protein in rabbit and affinity purified with Protein G beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	Biotin
-----------	--------

### Applications

---

Rabbit Anti-Alexa Fluor 555 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with biotin. This configuration enables flexible signal detection using streptavidin- or avidin-based systems in various assay formats.

The antibody may cross-react with related fluorophores such as iFluor® 555, ATTO 550, and Cy3. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows.

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