

## Rabbit Anti-Alexa Fluor 555 Antibody

Catalog Number: 46530

Unit Size: 1 mg

### 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, 0.05% Sodium Azide

### Antibody Properties

---

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

### Biological Properties

---

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Alexa Fluor 555 coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Applications

---

Rabbit Anti-Alexa Fluor 555 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated biomolecules. It offers high-affinity detection of Alexa Fluor® 555-labeled proteins, nucleic acids, or other tagged targets in immunoassays and imaging workflows.

The antibody may cross-react with spectrally similar fluorophores such as iFluor® 555, ATTO 550, and Cy3. It is compatible with standard secondary antibody systems and is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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