

## Rabbit Anti-Alexa Fluor 750 Antibody

Catalog Number: 46590

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 750 coupled to carrier protein

### Biological Properties

---

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Alexa Fluor 750 coupled carrier protein in rabbit and affinity purified with Protein G beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Applications

---

Rabbit Anti-Alexa Fluor 750 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules. It is ideal for detecting Alexa Fluor® 750-labeled antibodies, proteins, or probes in fluorescence-based applications.

The antibody may cross-react with structurally related fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. It is compatible with techniques including Western blotting, immunofluorescence microscopy, ELISA, and flow cytometry.

*Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*