

## Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 750-Labeled

Catalog Number: 46512

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	Alexa Fluor 488 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 488 coupled carrier protein in rabbit and affinity purified with Protein G beads. The antibody is conjugated with iFluor® 750 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

### Applications

---

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima around 749/775 nm. The direct labeling enables single-step detection of Alexa Fluor® 488-labeled molecules without the need for a secondary antibody.

The antibody may cross-react with structurally related green fluorophores such as iFluor® 488, ATTO 488, and rhodamine 110. With its long-wavelength emission, iFluor® 750 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.*