

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

Catalog Number: 46595

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, BSA

### Unit Details

---

Reconstitution Volume	0.1 mL ddH2O
-----------------------	--------------

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 750 coupled to carrier protein

### Biological Properties

---

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

### Spectral Properties

---

Conjugate	iFluor® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

### Applications

---

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 488, a bright green-emitting fluorophore with excitation/emission maxima of ~491/516 nm. This direct labeling format enables single-step detection of Alexa Fluor® 750 targets without secondary antibodies. The antibody may cross-react with structurally similar dyes such as iFluor® 750, ATTO 740, and DyLight® 755. Its spectral separation from Alexa Fluor® 750 makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*