

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

Catalog Number: 46513

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 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® 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 488 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 488, a green-emitting fluorophore with excitation around 490 nm. This direct labeling format allows sensitive, single-step fluorescence detection, minimizing background and improving workflow efficiency.

The antibody may cross-react with structurally similar dyes such as iFluor® 488, ATTO 488, and rhodamine 110. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.

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