

## Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 555-Labeled

Catalog Number: 46505

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 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 A beads. The antibody is conjugated with iFluor® 555 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

### Applications

---

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor<sup>®</sup> 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor<sup>®</sup> 488-conjugated targets and is directly labeled with iFluor<sup>®</sup> 555, a bright, orange-red fluorophore excited at 550–560 nm. This direct labeling allows for single-step detection of Alexa Fluor 488-tagged biomolecules without requiring a secondary antibody. The antibody may cross-react with similar green fluorophores such as iFluor<sup>®</sup> 488, ATTO 488, and rhodamine 110. Its emission in the orange-red spectrum enables spectral separation from Alexa Fluor<sup>®</sup> 488, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor<sup>®</sup> is a registered trademark of Thermo Fisher Scientific.*