

## Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 647-Labeled

Catalog Number: 46508

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 ddH <sub>2</sub> O
-----------------------	---------------------------

### 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® 647 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 647
Excitation Wavelength	656 nm
Emission Wavelength	670 nm

### Applications

---

Rabbit Anti-Alexa Fluor® 488 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. The direct labeling format allows for one-step detection of Alexa Fluor® 488-labeled molecules without the need for a secondary antibody. The antibody may cross-react with similar green fluorophores such as iFluor® 488, ATTO 488, and rhodamine 110. Its far-red emission makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.*