

# Rabbit Anti-Alexa Fluor 488 Antibody, Biotinylated

Catalog Number: 46520

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

## Spectral Properties

---

Conjugate	Biotin
-----------	--------

## Applications

---

Rabbit Anti-Alexa Fluor 488 Antibody, Biotinylated is a polyclonal antibody that specifically recognizes Alexa Fluor® 488-conjugated molecules and is directly conjugated to biotin. The biotin tag enables flexible signal amplification using streptavidin, or other avidin-based systems, making it suitable for diverse assay formats.

The antibody may cross-react with structurally related dyes such as iFluor® 488, ATTO 488, and rhodamine 110. Its broad binding and versatile detection makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence

detection.

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