

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, 0.2% (w/v) BSA

Unit Details

Reconstitution	0.1 mL ddH ₂ 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 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.