

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 350-Labeled

Catalog Number: 46502

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 ddH2O
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® 350 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 350, a bright, UV-excitable blue fluorophore. This format allows for direct detection of Alexa Fluor 488-labeled biomolecules without the need for a secondary antibody.

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

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