

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

Catalog Number: 46535

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 555 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 555 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 488 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 488, a bright green-emitting fluorophore with excitation/emission maxima around 491/517 nm. This direct labeling format eliminates the need for a secondary antibody, enabling fast and simplified detection of Alexa Fluor® 555-labeled biomolecules.

The antibody may cross-react with structurally related dyes such as iFluor® 555, ATTO 550, and Cy3. Its green emission makes it compatible with multicolor fluorescence assays that require distinct spectral separation from red and far-red fluorophores. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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