

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled

Catalog Number: 46542

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 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 G beads. The antibody is conjugated with iFluor® 750 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

Applications

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima of ~752/778 nm. This direct labeling enables single-step fluorescence detection without the need for secondary antibodies.

The antibody may cross-react with similar fluorophores such as iFluor® 555, ATTO 550, and Cy3. The near-infrared emission of iFluor® 750 allows for low-background detection and is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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