

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 555-Labeled

Catalog Number: 46628

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 350 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 350 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 555 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 555, a bright orange fluorophore with excitation/emission at around 555/565 nm. This direct labeling format enables one-step fluorescence detection without the need for a secondary antibody. The antibody may cross-react with related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its orange emission is spectrally distinct from Alexa Fluor® 350, making it ideal for multicolor imaging, flow cytometry, and other fluorescence-based assays. Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.