

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

Catalog Number: 46543

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 ddH ₂ O
-----------------------	---------------------------

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® 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 555 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation around 555 nm. This direct labeling format enables single-step detection, improving signal sensitivity and reducing background interference. The antibody may cross-react with structurally similar dyes such as iFluor® 555, ATTO 550, and Cy3. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation, and performing multiplex fluorescence assays. *Alexa Fluor® and Cy3 are registered trademarks of Thermo Fisher Scientific.*