

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

Catalog Number: 46633

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 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 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 350 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 350, a UV-excitabile blue fluorophore with excitation around 350 nm and emission near 442 nm. This direct labeling enables one-step fluorescence detection, reducing assay complexity and minimizing background signal.

The antibody may cross-react with structurally similar dyes such as iFluor® 350, AMCA, and Cascade Blue®. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in microscopy, ELISA, and flow cytometry.

Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.