

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 750-Labeled

Catalog Number: 46632

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 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® 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 350 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission at around 752/778 nm. This direct format allows single-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with structurally related fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its NIR emission provides high signal-to-noise ratios and is well-suited for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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