

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

Catalog Number: 46603

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 750 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 750 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 750 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation around 752 nm and emission near 778 nm. This direct labeling format allows single-step fluorescence detection with high sensitivity and minimal background.

The antibody may cross-react with structurally similar dyes such as iFluor® 750, ATTO 740, and DyLight® 755. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.

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