

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

Catalog Number: 46592

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 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® 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 750 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 350, a UV-excitabile blue fluorophore with excitation/emission of ~346/442 nm. This direct labeling format enables one-step detection of Alexa Fluor® 750-labeled targets without the need for secondary antibodies.

The antibody may cross-react with structurally related dyes such as iFluor® 750, ATTO 740, and DyLight® 755. Its blue emission is well separated from near-infrared signals, making it ideal for multiplex assays and low-background imaging workflows.

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