

Rabbit Anti-Cy7 Antibody, iFluor® 750-Labeled

Catalog Number: 46951

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	Cy7 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	Antibody is produced with immunogen in rabbit, affinity purified and 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-Cy7 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Cy7-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission around 752/778 nm. This direct labeling enables sensitive, single-step detection of Cy7-labeled molecules without secondary antibodies.

The antibody may cross-react with similar dyes such as Alexa Fluor® 750, DyLight® 755, and other far-red fluorophores. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in deep-tissue imaging, flow cytometry, and high-content analysis.

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