

Rabbit Anti-Cy7 Antibody, iFluor® 488-Labeled

Catalog Number: 46945

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 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	Liquid
Preparation	It is produced with Cy7 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 488 under optimal conditions.
Soluble In	Water
Application	ELISA, IF, FACS

Spectral Properties

Conjugate	iFluor 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-Cy7 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically recognizes Cy7-labeled targets and is directly conjugated to iFluor® 488, a bright green fluorophore with excitation/emission maxima of 491/516 nm. This direct labeling format enables single-step fluorescence detection and streamlined assay workflows without the need for secondary antibodies. The antibody may show cross-reactivity with other structurally related cyanine dyes. It is suitable for immunofluorescence microscopy, flow cytometry, and other fluorescence-based applications where spectrally separated detection of Cy7-labeled probes is required. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.