

Rabbit Anti-Cy3 Antibody, iFluor® 488-Labeled

Catalog Number: 46885

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	Cy3 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Cy3 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-Cy3 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Cy3-conjugated molecules and is directly labeled with iFluor® 488, a green-fluorescent dye with excitation/emission maxima of ~491/516 nm. This direct labeling format enables one-step detection of Cy3-labeled biomolecules without requiring a secondary antibody. The antibody may exhibit cross-reactivity with related cyanine based dyes. It is ideal for signal amplification, dye conjugation validation, and multicolor fluorescence detection in microscopy, flow cytometry, and other imaging applications.