

Rabbit Anti-Cy5 Antibody, iFluor® 555-Labeled

Catalog Number: 46918

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	Cy5 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Cy5 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 555 under optimal conditions.
Soluble In	Water
Application	ELISA, IF, FACS

Spectral Properties

Conjugate	iFluor 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-Cy5 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Cy5-conjugated molecules and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission maxima of ~557/570 nm. This direct labeling format supports one-step detection of Cy5-labeled biomolecules without the need for a secondary antibody. The antibody may exhibit cross-reactivity with related cyanine-based dyes. It is recommended for dye conjugation validation, signal amplification in low-abundance targets, and multicolor fluorescence applications in microscopy and flow cytometry.