

Rabbit Anti-Cy3 Antibody, iFluor® 647-Labeled

Catalog Number: 46888

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® 647 under optimal conditions.
Soluble In	Water
Application	ELISA, IF, FACS

Spectral Properties

Conjugate	iFluor 647
Excitation Wavelength	656 nm
Emission Wavelength	670 nm

Applications

Rabbit Anti-Cy3 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Cy3-conjugated molecules and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. This direct labeling format supports one-step detection of Cy3-labeled biomolecules without the need for a secondary antibody. The antibody may exhibit cross-reactivity with related cyanine based dyes. It is recommended for signal amplification, conjugation validation, and multicolor fluorescence applications in microscopy and flow cytometry.