

Rabbit Anti-Cy5 Antibody, iFluor® 647-Labeled

Catalog Number: 46921

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	Cy5 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	Antibody is produced with immunogen in rabbit, affinity purified and conjugated with iFluor® 647 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 647
Excitation Wavelength	656 nm
Emission Wavelength	670 nm

Applications

Rabbit Anti-Cy5 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Cy5-conjugated molecules and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission around 656/670 nm. This direct labeling format provides efficient, single-step fluorescence detection without the need for secondary antibodies.

The antibody may cross-react with similar dyes such as Alexa Fluor® 647, DyLight® 650, and ATTO 647. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays.

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