

Rabbit Anti-Cy5 Antibody

Catalog Number: 46910

Unit Size: 1 mg

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, 0.05% Sodium Azide

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Cy5 coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	Antibody is produced with immunogen in rabbit and affinity purified.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Rabbit Anti-Cy5 Antibody is a polyclonal antibody that specifically binds to Cy5-labeled molecules. It is designed to detect Cy5-conjugated proteins, oligonucleotides, or other biomolecules in fluorescence-based assays with high sensitivity.

The antibody may cross-react with structurally similar dyes including iFluor® 647, Alexa Fluor® 647, DyLight® 650, and ATTO 647. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and supporting multiplex fluorescence workflows in microscopy, flow cytometry, and blotting platforms.

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