

Rabbit Anti-Cy3 Antibody

Catalog Number: 46880

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	Cy3 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-Cy3 Antibody is a polyclonal antibody that specifically binds to Cy3-conjugated biomolecules. It is designed to recognize the cyanine dye Cy3 commonly used for fluorescent labeling in microscopy, flow cytometry, and molecular biology assays.

The antibody may also cross-react with structurally similar fluorophores such as Alexa Fluor® 555, ATTO 550, and iFluor® 555. It is ideal for signal amplification in low-abundance target detection, verifying Cy3 conjugation efficiency, and enabling multiplex fluorescence analysis in various platforms.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.