

Rabbit Anti-Cy7 Antibody

Catalog Number: 46940

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	Cy7 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-Cy7 Antibody is a polyclonal antibody that specifically binds to molecules labeled with the near-infrared dye Cy7. It is designed for detecting Cy7-conjugated antibodies, proteins, or nucleic acids in fluorescence-based applications.

The antibody may cross-react with spectrally similar fluorophores such as Alexa Fluor® 750, DyLight® 755, and iFluor® 750. It is ideal for signal amplification in low-abundance target detection, verifying conjugation efficiency, and enabling multiplexed fluorescence assays.

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