

Rabbit Anti-Cy5 Antibody, Biotinylated

Catalog Number: 46930

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 ddH ₂ O
-----------------------	---------------------------

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 Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Rabbit Anti-Cy5 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Cy5-conjugated molecules and is directly labeled with biotin. This format enables flexible signal detection using streptavidin- or avidin-based reporters in both colorimetric and fluorescence-based assays.

The antibody may cross-react with structurally related 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 supporting multiplexed analysis across ELISA, blotting, and imaging workflows.

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