

Rabbit Anti-Cy7 Antibody, Biotinylated

Catalog Number: 46960

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	Cy7 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-Cy7 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Cy7-conjugated targets and is directly labeled with biotin. This biotinylated format enables flexible signal detection using streptavidin- or avidin-based systems across a range of assay platforms.

The antibody may cross-react with other near-infrared dyes 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 performing multiplex fluorescence assays.

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