

Rabbit Anti-Texas Red Antibody, Biotinylated

Catalog Number: 46710

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, BSA

Unit Details

Reconstitution Volume	0.1 mL ddH2O
-----------------------	--------------

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Texas Red coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with ROX coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	Biotinylated
-----------	--------------

Applications

Rabbit Anti-Texas Red Antibody, Biotinylated is a polyclonal antibody that specifically binds to Texas Red®-conjugated molecules and is directly labeled with biotin. This configuration enables flexible detection using streptavidin- or avidin-linked systems, making it suitable for a variety of assay platforms. The antibody may cross-react with structurally similar fluorophores such as Alexa Fluor® 594 and ATTO 594. It is ideal for signal amplification in low-abundance target detection, verifying dye

conjugation efficiency, and performing multiplex fluorescence assays in applications such as microscopy, ELISA, and blotting. *Texas Red® and Alexa Fluor® are registered trademarks of Thermo Fisher Scientific.*