

## 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, 0.2% (w/v) BSA

### Unit Details

---

Reconstitution Volume 0.1 mL ddH<sub>2</sub>O

### 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 G beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### 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.*