

## Rabbit Anti-Texas Red Antibody, iFluor® 594-Labeled

Catalog Number: 46693

Unit Size: 100 ug

### Product Details

---

Storage Conditions 2-8°C with minimized light exposure. Do not freeze.

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 Texas Red coupled carrier protein in rabbit and affinity purified with Protein G beads. The antibody is conjugated with iFluor® 594 under optimal conditions.

Soluble In Water

Application IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate iFluor® 594

Excitation Wavelength 587 nm

Emission Wavelength 603 nm

### Applications

---

Rabbit Anti-Texas Red Antibody, iFluor® 594-Labeled is a polyclonal antibody that specifically binds to Texas Red®-conjugated molecules and is directly labeled with iFluor® 594, a red-emitting fluorophore with excitation around 590 nm. This format

enables single-step fluorescence detection and signal amplification, eliminating the need for secondary antibodies and reducing background in complex workflows.

The antibody may cross-react with similar fluorophores such as iFluor® 594 and Alexa Fluor® 594. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in applications like microscopy, flow cytometry, and ELISA.

*Texas Red® and Alexa Fluor® are registered trademarks of Thermo Fisher Scientific.*