

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. |
| 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 Texas Red coupled carrier protein in rabbit and affinity purified with Protein A 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.*