

## Rabbit Anti-Texas Red Antibody, Biotinylated

Catalog Number: 46710

Unit Size: 100 ug

### Product Details

---

|                    |                                            |
|--------------------|--------------------------------------------|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
|--------------------|--------------------------------------------|

|                 |                          |
|-----------------|--------------------------|
| Expiration Date | 12 months upon receiving |
|-----------------|--------------------------|

### Chemical Properties

---

|            |        |
|------------|--------|
| Appearance | Liquid |
|------------|--------|

|                  |     |
|------------------|-----|
| Molecular Weight | N/A |
|------------------|-----|

|            |     |
|------------|-----|
| Soluble In | PBS |
|------------|-----|

### Spectral Properties

---

|                       |     |
|-----------------------|-----|
| Excitation Wavelength | N/A |
|-----------------------|-----|

|                     |     |
|---------------------|-----|
| Emission Wavelength | N/A |
|---------------------|-----|

### 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.*