

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 555-Labeled

Catalog Number: 46658

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 | Alexa Fluor 594 coupled to carrier protein |

Biological Properties

| | |
|-------------|---|
| Stabilizer | BSA |
| Appearance | Solid |
| Preparation | It is produced with Alexa Fluor 594 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 555 under optimal conditions. |
| Soluble In | Water |
| Application | IF, ELISA, Flow Cytometry (FACS) |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor® 555 |
| Excitation Wavelength | 557 nm |
| Emission Wavelength | 570 nm |

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission at around 556/570 nm. This direct format allows single-step fluorescence detection without the need for a secondary antibody. The antibody may cross-react with structurally related dyes such as iFluor® 594, ATTO 594, and Texas Red®. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.
Alexa Fluor® and Texas Red® are registered trademarks of Thermo Fisher Scientific.