

## Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled

Catalog Number: 46542

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 | 757 nm |
|-----------------------|--------|

|                     |        |
|---------------------|--------|
| Emission Wavelength | 779 nm |
|---------------------|--------|

### Applications

---

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima of ~752/778 nm. This direct labeling enables single-step fluorescence detection without the need for secondary antibodies.

The antibody may cross-react with similar fluorophores such as iFluor® 555, ATTO® 550, and Cy3. The near-infrared emission of iFluor® 750 allows for low-background detection and is ideal for deep-tissue imaging, multicolor assays, and high-sensitivity applications.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.