

## Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 488-Labeled

Catalog Number: 46655

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® 488 under optimal conditions. |
| Soluble In  | Water   |
| Application | IF, ELISA, Flow Cytometry (FACS)  |

### Spectral Properties

---

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor® 488 |
| Excitation Wavelength | 491 nm      |
| Emission Wavelength   | 516 nm      |

### Applications

---

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated molecules and is directly labeled with iFluor® 488, a green-emitting fluorophore excited at around 490 nm. This direct format enables one-step fluorescence detection without requiring a secondary antibody. The antibody may cross-react with structurally similar dyes such as iFluor® 594, ATTO 594, and Texas Red®. Its green fluorescence is spectrally separated from orange-red dyes, making it ideal for multicolor assays including microscopy, ELISA, and flow cytometry. *Alexa Fluor® and Texas Red® are registered trademarks of Thermo Fisher Scientific.*