

## Rabbit Anti-Fluorescein [FITC] Antibody, Biotinylated

Catalog Number: 46770

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	FITC coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Coumarin coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	Biotinylated
-----------	--------------

### Applications

---

Rabbit Anti-Fluorescein [FITC] Antibody, Biotinylated is a polyclonal antibody that specifically binds to fluorescein-based dyes such as FITC and is directly labeled with biotin. This configuration enables flexible detection using streptavidin- or avidin-based systems in a variety of assay formats. The antibody may cross-react with structurally related dyes such as Alexa Fluor® 488, DyLight® 488, and other fluorescein analogs. It is ideal for signal amplification in detecting low-abundance FITC conjugates,

validating conjugation efficiency, and supporting multiplexed detection workflows. *Alexa Fluor®* and *DyLight®* are registered trademarks of Thermo Fisher Scientific.