

# Rabbit Anti-Fluorescein [FITC] Antibody, iFluor® 488-Labeled

Catalog Number: 46661

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 FITC coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with Biotin 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-Fluorescein [FITC] Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to fluorescein-based dyes such as FITC and is directly labeled with iFluor® 488, a green-emitting fluorophore with excitation around 490 nm. This labeling format allows for sensitive, one-step fluorescence detection of FITC-conjugated targets without requiring a secondary antibody. The antibody may cross-react with related dyes including Alexa Fluor® 488, DyLight® 488, and other xanthene-based fluorophores. It is ideal for signal amplification in detecting low-abundance targets, verifying fluorescein conjugation, and enabling multiplex fluorescence assays in applications such as flow cytometry, microscopy, and ELISA. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*