

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, 0.2% (w/v) BSA

Unit Details

Reconstitution Volume	0.1 mL ddH ₂ O
-----------------------	---------------------------

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 G 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.