

Rabbit Anti-Fluorescein [FITC] Antibody

Catalog Number: 46750

Unit Size: 1 mg

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, 0.05% Sodium Azide

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	FITC coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with FITC 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)

Applications

Rabbit Anti-Fluorescein [FITC] Antibody is a polyclonal antibody that specifically binds to fluorescein-based dyes such as FITC (fluorescein isothiocyanate), a widely used green fluorophore in molecular biology. This antibody enables detection of FITC-labeled biomolecules including antibodies, oligonucleotides, and proteins. It may cross-react with structurally related dyes like Alexa Fluor® 488, and DyLight® 488. The antibody is ideal for signal amplification in applications where FITC-labeled targets are present at low levels, and is suitable for use in various techniques. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*