

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

Catalog Number: 46513

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 488 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 488 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 488 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 488, a green-emitting fluorophore with excitation around 490 nm. This direct labeling format allows sensitive, single-step fluorescence detection, minimizing background and improving workflow efficiency. The antibody may cross-react with structurally similar dyes such as iFluor® 488, ATTO 488, and rhodamine 110. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications. *Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.*