

Rabbit Anti-Alexa Fluor 488 Antibody

Catalog Number: 46500

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	Alexa Fluor 488 coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Alexa Fluor 488 coupled carrier protein in rabbit and affinity purified with Protein G beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Rabbit Anti-Alexa Fluor 488 Antibody is a polyclonal antibody that binds to conjugates labeled with Alexa Fluor® 488 dye. It offers high-affinity binding, making it ideal for applications requiring signal amplification or dye conjugation verification.

The antibody may cross-react with similar fluorophores such as iFluor® 488, ATTO 488, and rhodamine 110. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.