

Rabbit Anti-Alexa Fluor 350 Antibody

Catalog Number: 46620

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 350 coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Alexa Fluor 350 coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Rabbit Anti-Alexa Fluor 350 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules. It is designed for detecting Alexa Fluor® 350-labeled antibodies, proteins, or other probes in fluorescence-based assays. The antibody may cross-react with related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.*