

Rabbit Anti-Alexa Fluor 647 Antibody

Catalog Number: 46560

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 647 coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Alexa Fluor 647 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 647 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules. It is ideal for detecting Alexa Fluor® 647-labeled antibodies, proteins, or probes. The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO 647 and DyLight® 650. It is ideal for applications requiring signal amplification or dye conjugation verification. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*