

Rabbit Anti-Alexa Fluor 647 Antibody, Biotinylated

Catalog Number: 46580

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, 0.2% (w/v) BSA

Unit Details

Reconstitution	0.1 mL ddH ₂ O
Volume	

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 647 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 647 coupled carrier protein in rabbit and affinity purified with Protein G beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	Biotin
-----------	--------

Applications

Rabbit Anti-Alexa Fluor 647 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with biotin. This configuration enables flexible signal detection using streptavidin- or avidin-based systems in a wide range of assay formats.

The antibody may cross-react with structurally similar fluorophores such as iFluor® 647, Cy5, ATTO 647, and DyLight® 650. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.