

Rabbit Anti-Alexa Fluor 594 Antibody, Biotinylated

Catalog Number: 46670

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 594 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 594 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 594 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated targets and is directly labeled with biotin. This configuration enables flexible signal detection using streptavidin- or avidin-linked systems in a wide range of assay formats.

The antibody may cross-react with structurally similar fluorophores such as iFluor® 594, ATTO 594, and Texas Red®. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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