

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 594-Labeled

Catalog Number: 46663

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, BSA

Unit Details

Reconstitution Volume	0.1 mL ddH2O
-----------------------	--------------

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 A beads. The antibody is conjugated with iFluor® 594 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 594
Excitation Wavelength	587 nm
Emission Wavelength	603 nm

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 594-Labeled is a polyclonal antibody that selectively targets Alexa Fluor® 594-conjugated biomolecules. It is directly conjugated to iFluor® 594, a bright red-emitting fluorophore with excitation/emission maxima of ~590/617 nm, enabling robust one-step detection of Alexa Fluor® 594-labeled targets in immunofluorescence, microscopy, and flow cytometry. This antibody may cross-react with spectrally similar dyes such as DyLight® 594, and ATTO 594. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*