

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 555-Labeled

Catalog Number: 46658

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 iFluor® 555 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission at around 556/570 nm. This direct format allows single-step fluorescence detection without the need for a secondary antibody.

The antibody may cross-react with structurally related dyes 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.