

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 488-Labeled

Catalog Number: 46655

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

Spectral Properties

Conjugate	iFluor® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated molecules and is directly labeled with iFluor® 488, a green-emitting fluorophore excited at around 490 nm. This direct format enables one-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with structurally similar dyes such as iFluor® 594, ATTO 594, and Texas Red®. Its green fluorescence is spectrally separated from orange-red dyes, making it ideal for multicolor assays including microscopy, ELISA, and flow cytometry.

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