

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 555-Labeled

Catalog Number: 46568

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 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 A 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 647 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with iFluor® 555, a bright orange-red fluorophore with excitation/emission of ~555/565 nm. This direct format enables one-step detection of Alexa Fluor® 647-labeled targets without the need for secondary antibodies. The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO 647, and DyLight® 650. The orange-red fluorescence of iFluor® 555 is spectrally distinct from far-red dyes, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*