

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 488-Labeled

Catalog Number: 46565

Unit Size: 100 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	PBS

Spectral Properties

Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is labeled with iFluor® 488, a green fluorescent dye with excitation/emission of ~491/516 nm. This direct labeling format allows 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. Its green fluorescence is spectrally distinct from far-red targets, enabling clean multiplex detection in immunofluorescence, ELISA, and flow cytometry workflows.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.