

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 350-Labeled

Catalog Number: 46562

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	345 nm
Emission Wavelength	450 nm

Applications

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is labeled with iFluor® 350, a bright, UV-excitable fluorophore with excitation/emission maxima of ~350/442 nm. This direct labeling format enables one-step fluorescence detection 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 spectrally distinct blue emission makes it ideal for multiplex applications where separation from far-red fluorophores is essential, supporting workflows in microscopy, ELISA, and flow cytometry.

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