

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 488-Labeled

Catalog Number: 46625

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 350 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 488, a bright green fluorophore with excitation/emission maxima of ~490/520 nm. This direct labeling allows for single-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its green fluorescence is well-separated from the blue emission of Alexa Fluor® 350, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.