

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

Catalog Number: 46502

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 488 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 350, a bright, UV-excitable blue fluorophore. This format allows for direct detection of Alexa Fluor 488-labeled biomolecules without the need for a secondary antibody.

The antibody may cross-react with structurally related dyes such as iFluor® 488, ATTO 488, and rhodamine 110. Its spectral separation from Alexa Fluor 488 makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific. iFluor® is a registered trademark of AAT Bioquest, Inc.