

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 555-Labeled

Catalog Number: 46505

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 488 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 555, a bright, orange-red fluorophore excited at 550–560 nm. This direct labeling allows for single-step detection of Alexa Fluor 488-tagged biomolecules without requiring a secondary antibody.

The antibody may cross-react with similar green fluorophores such as iFluor® 488, ATTO® 488, and rhodamine 110. Its emission in the orange-red spectrum enables spectral separation from Alexa Fluor® 488, making it ideal for multiplex fluorescence applications including flow cytometry, microscopy, and high-throughput screening.

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