

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 750-Labeled

Catalog Number: 46512

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	757 nm
Emission Wavelength	779 nm

Applications

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima around 749/775 nm. The direct labeling enables single-step detection of Alexa Fluor® 488-labeled molecules without the need for a secondary antibody.

The antibody may cross-react with structurally related green fluorophores such as iFluor® 488, ATTO® 488, and rhodamine 110. With its long-wavelength emission, iFluor® 750 is ideal for multiplexing alongside visible spectrum dyes, making this reagent suitable for fluorescence imaging, flow cytometry, and high-throughput assays.

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