

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 647-Labeled

Catalog Number: 46508

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	656 nm
Emission Wavelength	670 nm

Applications

Rabbit Anti-Alexa Fluor® 488 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. The direct labeling format allows for one-step detection of Alexa Fluor® 488-labeled molecules without the need for a secondary antibody.

The antibody may cross-react with similar green fluorophores such as iFluor® 488, ATTO® 488, and rhodamine 110. Its far-red emission makes it ideal for multiplex fluorescence applications by avoiding spectral overlap with green dyes, supporting workflows in immunofluorescence microscopy, 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.