

Rabbit Anti-Alexa Fluor 488 Antibody

Catalog Number: 46500

Unit Size: 1 mg

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	N/A
Emission Wavelength	N/A

Applications

Rabbit Anti-Alexa Fluor 488 Antibody is a polyclonal antibody that binds to conjugates labeled with Alexa Fluor® 488 dye. It offers high-affinity binding, making it ideal for applications requiring signal amplification or dye conjugation verification.

The antibody may cross-react with similar fluorophores such as iFluor® 488, ATTO® 488, and rhodamine 110. Its compatibility with techniques like immunofluorescence, Western blotting, ELISA, and flow cytometry makes it a versatile tool for imaging and analytical workflows.

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