

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

Catalog Number: 46513

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® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 488, a green-emitting fluorophore with excitation around 490 nm. This direct labeling format allows sensitive, single-step fluorescence detection, minimizing background and improving workflow efficiency.

The antibody may cross-react with structurally similar dyes such as iFluor® 488, ATTO® 488, and rhodamine 110. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.

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