

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 488-Labeled

Catalog Number: 46655

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 594 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated molecules and is directly labeled with iFluor® 488, a green-emitting fluorophore excited at around 490 nm. This direct format enables one-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with structurally similar dyes such as iFluor® 594, ATTO® 594, and Texas Red®. Its green fluorescence is spectrally separated from orange-red dyes, making it ideal for multicolor assays including microscopy, ELISA, and flow cytometry.

Alexa Fluor® and Texas Red® are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.