

Rabbit Anti-Texas Red Antibody, iFluor® 594-Labeled

Catalog Number: 46693

Unit Size: 100 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	
-----------------	--

Chemical Properties

Appearance	Liquid
------------	--------

Molecular Weight	N/A
------------------	-----

Soluble In	PBS
------------	-----

Spectral Properties

Excitation Wavelength	587 nm
-----------------------	--------

Emission Wavelength	603 nm
---------------------	--------

Applications

Rabbit Anti-Texas Red Antibody, iFluor® 594-Labeled is a polyclonal antibody that specifically binds to Texas Red®-conjugated molecules and is directly labeled with iFluor® 594, a red-emitting fluorophore with excitation around 590 nm. This format enables single-step fluorescence detection and signal amplification, eliminating the need for secondary antibodies and reducing background in complex workflows.

The antibody may cross-react with similar fluorophores such as iFluor® 594 and Alexa Fluor® 594. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in applications like microscopy, flow cytometry, and ELISA.

Texas Red® and Alexa Fluor® are registered trademarks of Thermo Fisher Scientific.