

Rabbit Anti-Texas Red Antibody

Catalog Number: 46680

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-Texas Red Antibody is a polyclonal antibody that specifically binds to molecules conjugated with Texas Red®, a red-emitting fluorophore commonly used in fluorescence imaging and flow cytometry. This antibody offers strong affinity for Texas Red®-labeled targets and can be used in a variety of detection formats.

The antibody may cross-react with similar fluorophores such as Alexa Fluor® 594 and ATTO 594. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection in applications such as microscopy, ELISA, and blotting.

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