

Rabbit Anti-Texas Red Antibody

Catalog Number: 46680

Unit Size: 1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.05% Sodium Azide

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Texas Red coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Texas Red coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

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.*