

Rabbit Anti-ROX Antibody

Catalog Number: 46720

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	ROX 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. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Rabbit Anti-ROX Antibody is a polyclonal antibody that specifically binds to ROX, a red-emitting dye commonly used as a reference fluorophore in qPCR and probe-based assays. It enables sensitive detection and validation of ROX-conjugated oligonucleotides, probes, and labeled biomolecules. The antibody may cross-react with structurally similar rhodamine derivatives. It is ideal for verifying dye conjugation, enhancing detection sensitivity in low-abundance targets, and supporting multiplexed fluorescence applications in nucleic acid and protein-based assays.