

Rabbit Anti-ROX Antibody

Catalog Number: 46720

Unit Size: 1 mg

Product Details

Storage Conditions	Refrigerated (2-8 °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-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.