

Rabbit Anti-ROX Antibody, Biotinylated

Catalog Number: 46740

Unit Size: 100 ug

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-ROX Antibody, Biotinylated is a polyclonal antibody that specifically binds to ROX conjugated molecules and is directly labeled with biotin. This format enables flexible detection using streptavidin- or avidin-based signal amplification systems in non-PCR applications.

This is suitable for confirming ROX conjugation to probes or biomolecules, or for use in fluorescence-based detection assays such as microscopy, ELISA, and bead-based multiplex assays.