

## Rabbit Anti-ROX Antibody, Biotinylated

Catalog Number: 46740

Unit Size: 100 ug

### 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, BSA

### Unit Details

---

Reconstitution Volume	0.1 mL ddH <sub>2</sub> O
-----------------------	---------------------------

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	ROX coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with FITC coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	Biotinylated
-----------	--------------

### 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.