

---

## **cry3Aa Polyclonal Antibody, HRP Conjugated**

(Catalog # A54550)

---

### **Background**

Promotes colloid osmotic lysis by binding to the midgut epithelial cells of Coleoptera.

### **Description**

cry3Aa Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

### **Formulation**

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### **Specificity**

Bacillus thuringiensis tenebrionis

### **Isotype**

IgG

### **Uniprot ID**

P0A379

### **Purification**

>95%, Protein G purified

### **Immunogen**

Recombinant Bacillus thuringiensis subsp. Tenebrionis Pesticidal crystal protein cry3Aa protein (58-644AA)

### **Storage**

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### **Alternative Names**

73 kDa crystal protein Crystalline entomocidal protoxin Insecticidal delta-endotoxin CryIIIa(a) cry3Aa bt13, cry3A, cryC, cryIIIa, cryIIIa(a)

### **Application**

ELISA; Recommended dilution: ELISA 1:500-1:1000

---