

---

## Riboflavin-Binding Protein Polyclonal Antibody

(Catalog #A55691)

---

### Background

Required for the transport of riboflavin to the developing oocyte.

### Description

Riboflavin-Binding Protein Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

chicken

### Isotype

IgG

### Uniprot ID

P02752

### Purification

Protein G purified

### Immunogen

Recombinant Chicken Riboflavin-binding protein (18-225AA)

### 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

RBP,

### Application

ELISA; Recommended dilution: ELISA 1:1000-1:5000

---