

---

# Ovomucoid Polyclonal Antibody

(Catalog # A57565)

---

## Description

Ovomucoid 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

P01005

## Purification

>95%, Protein G purified

## Immunogen

Chicken Ovomucoid (Native Protein)

## Storage

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

## Application

ELISA

---