

---

## Tlr7 Polyclonal Antibody, Biotin Conjugated

(Catalog # A53660)

---

### Description

Tlr7 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

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

### Specificity

Mouse

### Isotype

IgG

### Uniprot ID

P58681

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Mouse Toll-like receptor 7 protein (27-348AA)

### 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

Tlr7

### Application

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

---