

---

## groL Polyclonal Antibody, HRP Conjugated

(Catalog # A50858)

---

### Background

Prevents misfolding and promotes the refolding and proper assembly of unfolded polypeptides generated under stress conditions.

### Description

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

### Formulation

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

### Specificity

Yersinia pestis

### Isotype

IgG

### Uniprot ID

A4TRR0

### Purification

Protein G purified

### Immunogen

Recombinant Yersinia pestis 60 kDa chaperonin protein (1-548AA)

### 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

GroEL protein;Protein Cpn60;groL;groEL

### Application

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

---