

## fabD Polyclonal Antibody

(Catalog # A50961)

---

### Description

fabD Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Staphylococcus aureus

### Isotype

IgG

### Uniprot ID

Q5HGK3

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Staphylococcus aureus Malonyl CoA-acyl carrier protein transacylase protein (1-308AA)

### 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

MCT, fabD, SACOL1244

### Application

ELISA

---