

---

## entC1 Polyclonal Antibody, HRP Conjugated

(Catalog #A50850)

---

### Background

Staphylococcal enterotoxins cause the intoxication staphylococcal food poisoning syndrome. The illness characterized by high fever, hypotension, diarrhea, shock, and in some cases death.

### Description

entC1 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

### Formulation

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

### Specificity

Staphylococcus aureus

### Isotype

IgG

### Uniprot ID

P01553

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Staphylococcus aureus Enterotoxin type C-1 protein (28-266AA)

### 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

SEC1, entC1

### Application

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

---