

---

## sspB Polyclonal Antibody, HRP Conjugated

(Catalog #A54566)

---

### Background

Cysteine protease able to degrade elastin, fibrogen, fibronectin and kininogen. Exhibits a strong preference for substrates where arginine is preceded by a hydrophobic amino acid. Promotes detachment of primary human keratinocytes. Along with other extracellular proteases is involved in colonization and infection of human tissues

### Description

sspB 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

Q99V46

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Staphylococcus aureus Staphopain B protein (220-393AA)

### 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

Staphylococcal cysteine proteinase B Staphylopain B sspB SAV1047

### Application

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

---