

---

## fim2 Polyclonal Antibody

(Catalog # A54740)

---

### Background

*Bordetella pertussis* is the causative agent of whooping cough. An essential step in the disease process is the attachment of the bacteria to the ciliated epithelium of the respiratory tract, enabling the organism to resist normal host-clearance mechanisms. It is unclear which bacterial cell surface component are responsible for adherence but the fimbriae of *B. pertussis* are prime candidates for being involved in this process.

### Description

fim2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

*Bordetella pertussis*

### Isotype

IgG

### Uniprot ID

P05788

### Purification

>95%, Protein G purified

### Immunogen

Recombinant *Bordetella pertussis* Serotype 2 fimbrial subunit protein (27-207AA)

### 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

fim2, BP1119

### Application

ELISA

---