

**B. pertussis Filamentous Hemagglutinin (FHA)**

Strain Tohama I  
native, Broth culture

| Cat. No.    | Amount |
|-------------|--------|
| PR-BA120F-S | 250 µg |
| PR-BA120F-L | 1 mg   |

**For general laboratory use.**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -80 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Form:** liquid

**Culture system:** Broth culture

**Protein concentration**

Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.

**Preservation**

50 % Glycerol

**Analysis**

ELISA and SDS-PAGE

**Preparation**

Pertussis FHA, highly purified from Bordetella pertussis culture supernatant. Pertussis FHA is presented in Tris buffer pH 8.0, 0.5 M NaCl and 50 % glycerol.

**Inactivation**

Whole B. pertussis organisms are removed by sterile filtration. Inactivity is confirmed by attempted growth under original culture conditions. Since no procedure can guarantee absolute sterility, the reagent should be handled with appropriate safety precautions.

**Agent**

Bordetella pertussis, Strain Tohama I