

---

## tetX Polyclonal Antibody

(Catalog # A50817)

---

### Background

Tetanus toxin acts by inhibiting neurotransmitter release. It binds to peripheral neuronal synapses, is internalized and moves by retrograde transport up the axon into the spinal cord where it can move between postsynaptic and presynaptic neurons. It inhibits neurotransmitter release by acting as a zinc endopeptidase that catalyzes the hydrolysis of the '76-Gln-/-Phe-77' bond of synaptobrevin-2.

### Description

tetX Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Clostridium tetani

### Isotype

IgG

### Uniprot ID

P04958

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Clostridium tetani Tetanus toxin protein (123-573AA)

### 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

Tentoxylisin, Tetanus toxin chain L, Tetanus toxin chain H, CTC\_p60, tetX

### Application

ELISA

---