
botB Polyclonal Antibody

(Catalog #A50736)

Background

Botulinum 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 cleaves the '76-Gln-/Phe-77' bond of synaptobrevin-2.

Description

botB Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Clostridium botulinum

Isotype

IgG

Uniprot ID

P10844

Purification

>95%, Protein G purified

Immunogen

Recombinant Clostridium botulinum Botulinum neurotoxin type B protein (1-427AA)

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

BoNT/B, Bontoxilysin-B, botB

Application

ELISA
