
ptxA Polyclonal Antibody

(Catalog #A53718)

Background

S1 is an NAD-dependent ADP-ribosyltransferase, which plays a crucial role in the pathogenesis of *B. pertussis* causing disruption of normal host cellular regulation. It catalyzes the ADP-ribosylation of a cysteine in the alpha subunit of host heterotrimeric G proteins. In the absence of G proteins it also catalyzes the cleavage of NAD⁺ into ADP-ribose and nicotinamide. It irreversibly uncouples the G-alpha GTP-binding proteins from their membrane receptors.

Description

ptxA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Bordetella pertussis

Isotype

IgG

Uniprot ID

P04977

Purification

Protein G purified

Immunogen

Recombinant *Bordetella pertussis* Pertussis toxin subunit 1 protein (35-269AA)

Storage

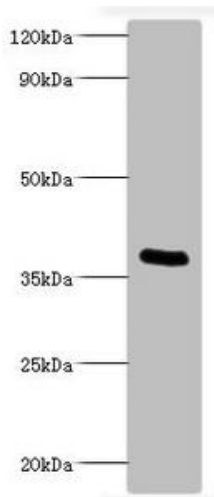
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Islet-activating protein S1 NAD-dependent ADP-ribosyltransferase ptxA BP3783

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: Bordella pertussis pertussis toxin subunit 1
antibody at 2ug/ml + recombinant Bordella pertussis
pertussis toxin subunit 1 100ng

Secondary

Goat polyclonal to rabbit IgG at 1/1000 dilution

Predicted band size: 36 kDa

Observed band size: 36 kDa