## Pertussis toxin (Bordetella pertussis)

Ultra-pure pertussis toxin delivering optimal antigenicity

Pertussis toxin consists of an enzymatically active A protomer subunit (S-1) which possesses both NAD<sup>+</sup> glycohydrolase and ADP-ribosyltransferase activities and a B oligomer subunit (S-2, S-3, S-4, and S-5) which is responsible for attachment of the native toxin to eukaryotic cell surfaces. Pertussis toxin uncouples G proteins from receptors by ADP ribosylating a cysteine near the carboxy-terminus of the A subunit.

Citations: 35

View Online »

**Ordering Information** 

Order Online »

BML-G100-0050

50µg

Manuals, SDS & CofA

View Online »

- Higher specificity and sensitivity
- Minimal lot-to-lot variation
- Reproducible scale-up for bulk quantities
- Lyophilized format for extended shelf-life

## **Handling & Storage**

**Use/Stability** 

As indicated on product label or CoA when stored as recommended. This preparation is lyophilized, aseptically packaged and sealed under vacuum. Do not vortex. Store at 4°C prior to and following reconstitution. Do not freeze. This toxin is considered to be biologically active. The lyophilized preparation will remain active for at least one year. After reconstitution with sterile water this product should remain active for at least six months if stored at 4°C. Each vial when reconstituted to 500 µl with sterile distilled water, contains 50.0 µg of pertussis toxin in 10 mM sodium phosphate buffer with 50 mM sodium chloride at pH 7.0. The toxin is not completely soluble and the resulting suspension should be made uniform by gentle mixing, not vortexing, prior to withdrawing aliquots. Do not sterile filter. For use with purified G proteins, the toxin must be structurally converted to artificially create a form of the toxin that interacts directly with G proteins. See Kaslow et al. (1987)This product should be used only by technically qualified personnel. Good laboratory technique should be employed: Avoid mouth pipetting; wear gloves and other protective clothing; avoid inhalation; avoid contact with skin or open wounds; thoroughly wash any area of the body that comes in contact with the toxin with soap and water; wash eyes with water for fifteen minutes; exercise care when handling in hypodermic needles to avoid inadvertent intravenous or subcutaneous injection.

**Handling** Aseptically packed, sealed under vacuum. Do not vortex. Do not freeze.

Long Term Storage +4°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Islet-activating protein, PTX, Holotoxin

**CAS** 70323-44-3

Formulation Lyophilized. Containing sodium phosphate and sodium

chloride.

Purity ≥95%

Purity Detail Purified by SDS-PAGE.

**Source** From *Bordetella pertussis*, strain 165.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com