

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⁺ 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. Activation Status: Please note that this product is supplied in a non-activated state. If your system requires activation, refer to Kaslow et al. (1987) Biochemistry 26, 123-127. 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. This product should be used only by technically qualified personnel. Good laboratory technique must be employed. This product should be used only by technically qualified personnel. Refer to the Safety Data Sheet (SDS) for complete handling, safety, and first aid information.
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 <i>Bordetella pertussis</i> , strain 165.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)