

Bordetella pertussis toxin

High purity toxin with minimal lot-to-lot variation

Native protein.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-G101-0050	50µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)

- High purity, native protein available as >98% pure
- Minimal lot-to-lot variation ensures reproducible results
- 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. 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. Resuspended toxin should be used within one month. 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.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Islet-activating protein, PTX
Appearance	White powder.
Application Notes	Antibody capture, ELISA, LINE assay, cytotoxicity studies
CAS	70323-44-3
Formulation	Lyophilized. Contains 0.05M sodium phosphate and 0.5M sodium chloride.
Purity	≥98%
Purity Detail	Purified protein.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.

From *Bordetella pertussis*, strain NCTC 10739.