

# C. difficile toxin B

Extremely potent cytotoxin.

Citations: 9

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-G150-0050	50µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)

- High potency retained following inactivation
- Minimal lot-to-lot variation
- Lyophilized for ease of handling and long term stability

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Clostridium difficile toxin B
Appearance	Clear residue.
Biological Activity	2 pg/ml by verocytotoxicity assay (lowest point at which cell rounding is visible)
Formulation	Lyophilized. Contains 0.05M HEPES, 0.15M sodium chloride and 5% sucrose.
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified protein.
Reconstitution	Reconstitute in 250µl sterile distilled water.
Source	From <i>Clostridium difficile</i> , strain VPI10463.
UniProt ID	P18177



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-eu@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)