

C. difficile toxin B from ribotype 027, (native)

Highly purified preparation from
hypervirulent ribotype 027

Potent enterotoxin.

Ordering Information

[Order Online »](#)

ENZ-PRT272-0002	2µg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

- High potency
- Minimal lot to lot variation
- Lypophilized for ease of handling and long term stability

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Clostridium difficile toxin B, tcdB, toxB
Formulation	Lyophilized. Contains 0.05M HEPES, 0.15M sodium chloride and 5% sucrose.
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified protein.
Reconstitution	Reconstitute in 7µl sterile distilled water.
Source	From <i>Clostridium difficile</i> , strain R20291 (Stoke Mandeville) ribotype 027.
UniProt ID	P18177

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-eu@enzolifesciences.com)

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