

C. difficile toxin A from ribotype 078, (native)

Highly purified preparation from
hypervirulent ribotype 078

Potent enterotoxin.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

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

Handling & Storage

Long Term Storage -80°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Clostridium difficile toxin A, tcdA, toxA
Formulation	Lyophilized. Contains 0.05M HEPES, 0.15M sodium chloride and 5% sucrose.
Purity	>90% (SDS-PAGE)
Purity Detail	Purified protein.
Reconstitution	Reconstitute in 17µl sterile distilled water.
Source	From <i>Clostridium difficile</i> , strain 210708 ribotype 078.
UniProt ID	P16154

Last modified: May 29, 2024



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-be@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