

Toxin A (Clostridium difficile)

Cytotoxic and enterotoxic

Has potent cytotoxic and enterotoxic properties. Translocates to the cytosol of target cells and glucosylates small GTP-binding proteins such as Rho, Rac, and Cdc42. Induces actin condensation, cell death and opening of the tight junctions. Triggers release of pro-inflammatory cytokines from intestinal epithelial cells probably via activation of MAP kinases.

Citations: 7

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 2 months at +4°C after reconstitution in sterile water at a concentration of 0.2 mg/ml.
Handling	Do not vortex. Do not freeze.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Activity	Each lot is tested to confirm binding activity to fresh rabbit red blood cells using a hemagglutination assay. The lowest concentration producing a positive reaction was ~50µg/ml.
Formulation	Lyophilized from 50mM TRIS, pH 7.5, containing 50mM sodium chloride and 0.1% trehalose.
MW	~270kDa
Purity Detail	When examined on 3-8% SDS-PAGE gels, this protein migrates as a single major band with an apparent molecular weight of 300,000 Da run. Densitometric analysis estimates the purity of the product as >80%.
Reconstitution	Reconstitute with sterile distilled water.
Source	Isolated from <i>Clostridium difficile</i> .
Technical Info / Product Notes	Protein concentration is determined by the Bradford method using BSA as a standard.
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