

# Toxin A (Clostridium difficile)

Cytotoxic and enterotoxigenic

Has potent cytotoxic and enterotoxigenic 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 »](#)

Ordering Information

[Order Online »](#)

ALX-630-100-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Do not vortex. Do not freeze.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Activity</b>	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.
<b>Formulation</b>	Lyophilized from 50mM TRIS, pH 7.5, containing 50mM sodium chloride and 0.1% trehalose.
<b>MW</b>	~270kDa
<b>Purity Detail</b>	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%.
<b>Reconstitution</b>	Reconstitute with sterile distilled water.
<b>Source</b>	Isolated from <i>Clostridium difficile</i> .
<b>Technical Info / Product Notes</b>	Protein concentration is determined by the Bradford method using BSA as a standard.
<b>UniProt ID</b>	P16154



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)