

Tetanus toxoid

Inactivation with formaldehyde leads to non-toxic form of tetanus toxin.

Produced from tetanus toxin by inactivation with formaldehyde. The resulting toxoid is essentially non-toxic.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-630-108-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 6 months when stored at -20°C.
Handling	Sealed under vacuum. Do not exceed one freeze/thaw cycle. Do not vortex. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Formaldehyde inactivated tetanus toxin
Concentration	0.5µg/µl after reconstitution.
Formulation	Lyophilized.
MW	~150kDa
Quality Control	Immunoprecipitation; concentration determination by UV absorbance.
Reconstitution	Reconstitute with 200µl distilled water to get a 0.5µg/µl solution in 100mM histidine, 100mM sodium chloride and 5% trehalose, pH 7.0.
UniProt ID	P04958

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)