

PRODUCT INFORMATION

Tetanus Toxin C-Fragment from *Clostridium tetani*

Item No. 23222

Overview and Properties

Contents:	Each vial, when reconstituted with 50 µl of sterile purified water, contains 10 µg of tetanus toxin C-fragment in 0.01 M sodium phosphate, pH 7.5
Storage:	4°C (as supplied)
Stability:	≥2 years
Protein	
Concentration:	Protein concentration was determined using a modified Bradford assay with ovalbumin as the standard

Special Conditions: Freezing is not recommended

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Tetanus toxin C-fragment from *C. tetani* is the C-terminal end of tetanus toxin that binds to gangliosides on presynaptic terminals at the neuromuscular junction, allowing tetanus toxin to be taken up into neurons.¹

Reference

1. Montecucco, C., Schiavo, F., Poulain, B., *et al.* Tetanus toxin is a zinc protein and its inhibition of neurotransmitter release and protease activity depend on zinc. *EMBO J.* **11(10)**, 3577-3583 (2016).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/09/2018

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM