

PRODUCT INFORMATION



Tetanus Toxin (1,301-1,315) (*C. tetani*)

Item No. 43402

CAS Registry No.: 158964-53-5

Formal Name: L-cysteinyl-L- α -aspartyl-L-tryptophyl-L-tyrosyl-L-phenylalanyl-L-valyl-L-prolyl-L-threonyl-L- α -aspartyl-L- α -glutamylglycyl-L-tryptophyl-L-threonyl-L-asparaginyl-L-aspartic acid

H—Cys—Asp—Trp—Tyr—Phe—Val—Pro—Thr—Asp—Glu—

Peptide Sequence: CDWYFVPTDEGWTND-OH

Gly—Trp—Thr—Asn—Asp—OH

MF: $C_{84}H_{106}N_{18}O_{28}S$

FW: 1,847.9

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

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

Laboratory Procedures

Tetanus toxin (1,301-1,315) (*C. tetani*) is supplied as a solid. A stock solution may be made by dissolving the tetanus toxin (1,301-1,315) (*C. tetani*) in the solvent of choice. Tetanus toxin (1,301-1,315) (*C. tetani*) is soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tetanus toxin (1,301-1,315) (*C. tetani*) is a C-terminal peptide fragment of tetanus toxin heavy chain from *C. tetani*.¹

Reference

1. Fischer, P.M. and Howden, M.E. Synthetic peptide antigens of tetanus toxin. *Mol. Immunol.* **31**(15), 1141-1148 (1994).

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, 07/21/2025

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