PRODUCT INFORMATION



Tetanus Toxoid

Item No. 19659

Overview and Properties

Source:	Each vial, when reconstituted with 50 μ l of water, contains 25 μ g of tetanus toxoid in 100 mM histidine, 100 mM NaCl, pH 7.0, and 5% trehalose
Storage:	2-8°C (as supplied); after reconstitution, aliquot, and freeze
Stability:	As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly
Protein	
Concentration:	Determined by absorbance at 280 nm using an extinction coefficient of 1.24 for a concentration of 1 mg/ml
Special conditions:	Handle gently; do not vortex

Description

Tetanus toxoid is prepared by formaldehyde inactivation of the neurotoxin, tetanus toxin. Formulations containing tetanus toxoid have been used as vaccines to produce immunity to tetanus, the disease caused by C. tetani infection.¹

Reference

1. Gupta, R.K. and Siber, G.R. Adjuvants for human vaccines--current status, problems and future prospects. Vaccine 13(14), 1263-1276 (1995).

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.

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