

PRODUCT INFORMATION



Diphtheria Toxoid from *Corynebacterium diphtheriae* Item No. 23219

Overview and Properties

Contents:	Each vial, when reconstituted to 1.0 ml with water, contains 500 µg of diphtheria toxoid in 0.01 M sodium phosphate, pH 7.5, containing 0.1 mg/ml sorbitol
Storage:	2-8°C (as supplied); after reconstitution store at -20°C
Stability:	≥4 years
Protein	
Concentration:	Determined by absorbance at 280 nm using an extinction coefficient of 1.18 for a concentration of 1 mg/ml

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

Description

Diphtheria toxoid is a formaldehyde inactivated form of diphtheria toxin that is non-specifically cross linked and subject to rearrangement.¹ Formulations containing diphtheria toxoid have been used as adjuvants in vaccines to produce immunity to diphtheria, the disease caused by *C. diphtheriae* infection.²

References

1. Jiang, W. and Schwendeman, S.P. Formaldehyde-mediated aggregation of protein antigens: Comparison of untreated and formalinized model antigens. *Biotechnol. Bioeng.* **70(5)**, 507-517 (2000).
2. Hansen, J., Zhang, L., Klein, N.P., et al. Post-licensure safety surveillance study of routine use of quadrivalent meningococcal diphtheria toxoid conjugate vaccine. *Vaccine* **35(49 Pt. B)**, 6879-6884 (2017).

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, 01/14/2026

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