

PRODUCT INFORMATION



Anti-Cholera Toxin B Subunit (goat)

Item No. 19655

Overview and Properties

Contents:	Each vial, when reconstituted with 0.1 ml of water, contains 0.1 ml of goat antiserum to cholera toxin B subunit (choleragenoid). Sodium azide (0.1%) has been added as a preservative
Form:	Solid
Storage:	4°C (as supplied); After reconstitution, aliquot, and freeze; Minimize freeze/thaw cycles
Stability:	≥2 years
Applications:	The greatest dilution of this lot, which forms an immunoprecipitin band against a 0.5 mg/ml solution of cholera toxin B subunit, is 1:4
Special Conditions:	Handle gently; do not vortex

Description

Anti-cholera toxin B subunit (goat) is a polyclonal anti-choleragenoid from goat that is suitable for use in toxin neutralization and binding assays or for immunohistochemical studies in conjunction with the cholera toxin B subunit. The pentameric B subunit of cholera toxin (Item No. 19654) binds with high efficiency to G_{M1} monosialoganglioside cell membrane receptors found on mammalian cell surfaces and facilitates entrance of the A subunit into cells.¹ The cholera toxin B subunit is often used for retrograde track tracing studies in neuroscience research.²

References

1. Cho, J.A., Chinnapen, D.J.F., Amar, E., *et al.* Insights on the trafficking and retro-translocation of glycosphingolipid-binding bacterial toxins. *Front. Cell. Infect. Microbiol.* **2(51)** (2012).
2. Baldauf, K.J., Royal, J.M., Hamorsky, K.T., *et al.* Cholera toxin B: One subunit with many pharmaceutical applications. *Toxins (Basel)* **7(3)**, 974-996 (2015).

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, 11/21/2022

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