



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Cholera Toxin Subunit B (Recombinant)

Recombinant cholera toxin subunit B (CTB) purified from *E. coli* as a single, non-glycosylated polypeptide chain. Free of stabilizers and ready to conjugate.



Product attributes

Probe cellular localization	Membrane/cell surface
Cell permeability	Membrane impermeant
Toxin	Cholera toxin
Storage Conditions	Store at -10 to -35 °C

Product Description

Biotium's recombinant cholera toxin subunit B (CTB) is purified from *E. coli* as a single, non-glycosylated polypeptide chain. Cholera toxin is the symptom-causing toxin produced by the bacteria *Vibrio cholerae* during cholera infection.

The active holo-toxin from Cholera is composed of a hexamer. Subunit A is the toxic enzymatic subunit present in one copy per toxin. Subunit B is the receptor binding subunit that is found as a pentamer in each toxin and is non-toxic, making it useful for cell biological studies. CTB is used as a macrophage stimulant (1) and a vaccine adjuvant (2). CTB also been used to determine endotoxin sensitivity of transgenic mice expressing human caspase-4 (3).

This purified recombinant CTB is free of stabilizers and ready to conjugate after reconstitution. Biotium also offers [CF® Dye labeled cholera toxin subunit B](#) that may be used as a lipid raft marker and endocytic tracer for imaging live or fixed cells.

References

1. Ther Adv Vaccines 3(5-6), 155-163 (2015), [DOI: 10.1177/2051013615613473](https://doi.org/10.1177/2051013615613473)
2. Infect Immun. 75(6), 3150-3159 (2007), [DOI: 10.1128/iai.00581-06](https://doi.org/10.1128/iai.00581-06)
3. J Immunol. 193(1), 335-343 (2014), [DOI: 10.4049/jimmunol.1303424](https://doi.org/10.4049/jimmunol.1303424)

This datasheet was generated on June 19, 2026 at 06:58:48 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cholera-toxin-subunit-b-recombinant/>