

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

Call us: 800-304-5357

Probe cellular localization	Membrane/cell surface
Cell permeability	Membrane impermeant
Toxin	Cholera toxin

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
 2. Infect Immun. 75(6), 3150-3159 (2007), DOI: 10.1128/iai.00581-06
 3. J Immunol. 193(1), 335-343 (2014), DOI: 10.4049/jimmunol.1303424

This datasheet was generated on November 29, 2025 at 08:23:06 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cholera-toxin-subunit-b-recombinant/