

Cholera toxin (Vibrio cholerae), (azide free)

G protein activator

May require a license for import, please [contact us](#) for more information.

Cholera toxin consists of a single A subunit surrounded by five B subunits. The B subunits are responsible for the attachment of the native toxin to ganglioside GM1 on mammalian cell surfaces, and facilitate translocation of the A subunit. The A subunit catalyzes the ADP-ribosylation of an arginine residue on the α subunit of G proteins, reducing intrinsic GTPase activity and activating the α subunit. ADP-ribosylation of other proteins such as human red cell Ca-ATPase has also been reported. Cholera toxin must interact with ADP-ribosylation factor (ARF) for maximal activity.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-G117-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Handle gently and do not vortex. Store at +4°C prior to and following reconstitution. Do not freeze.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

CAS	9012-63-9
Formulation	Lyophilized from 0.05M Tris, pH 7.5, containing 200 mM NaCl, and 1 mM Na ₂ EDTA.
Reconstitution	Reconstitute with 0.5ml sterile water to a 2mg/mL solution.
Source	From <i>Vibrio cholerae</i> .
UniProt ID	P01555 (A subunit), P01556 (B subunit)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)