C. difficile toxin B

Extremely potent cytotoxin.

Citations: 9

View Online »

Ordering Information

Order Online »

BML-G150-0050

50µg

Manuals, SDS & CofA

View Online »

- High potency retained following inactivation
- Minimal lot-to-lot variation
- Lyophilized for ease of handling and long term stability

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling After reconstitution, prepare aliquots and store at -80°C.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Clostridium difficile toxin B

Appearance Clear residue.

Biological Activity 2 pg/ml by verocytotoxicity assay (lowest point at which

cell rounding is visible)

Formulation Lyophilized. Contains 0.05M HEPES, 0.15M sodium

chloride and 5% sucrose.

Purity ≥95% (SDS-PAGE)

Purity Detail Purified protein.

Reconstitution Reconstitute in 250µl sterile distilled water.

Source From Clostridium difficile, strain VPI10463.

UniProt ID P18177

Last modified: June 13, 2024



eu@enzolifesciences.com