

Data Sheet

TGFβ1, Latent

Human, recombinant

Catalog #: 90901-1

Lot#: 141016

Formulation: Lyophilized from a 0.2 μM filtered, aqueous solution

Reconstitution: Supplied as a lyophilized solid. Reconstitute in sterile PBS with 0.1% BSA.

Stability: ≥12 months at -80°C (as a solid). Store reconstituted protein in aliquots at -20°C to -80°C for up to 6 months. Avoid freeze/thaw cycles.

References:

1. Derynck, R. *et al.* (1985). *Nature* **316(6030)**:701-5.
2. Bourdrel L., *et al.* (1993). *Protein Expr. Purif.* **4**:130-140.
3. Munger, J.S., *et al.* (1997). *Kidney Int.* **51**:1376-1382.

Description:

Human TGFβ1 (GenBank Accession No. NM_000660), a.a. 30-390 (end) (Cys33Ser), N-terminal His-tag, disulfide-linked homodimers of TGFβ1 LAP and mature TGFβ1, MW= ~75 & 25 kDa under non-reducing conditions.

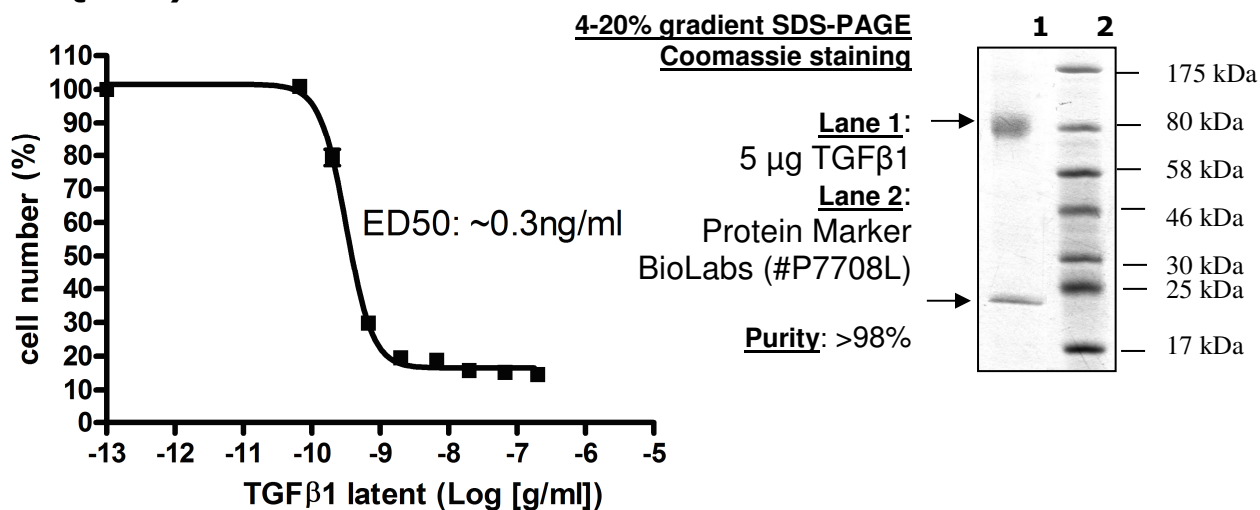
Source: FreeStyle 293-F cells.

Biological Activity: measured by TGFβ1's ability to inhibit mIL-4-dependent proliferation of HT-2 mouse helper T cells. The ED50 is <0.5 ng/ml.

Endotoxin Level: <1 EU/μg.

Application: Useful for cell culture and for the study of signaling and apoptotic pathways.

Quality Assurance



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com