

## Data Sheet

### TGFβ1, Active (Lyophilized)

Human, recombinant

**Catalog #:** 90900

**Lot#:** 180525

**Original concentration:**  
0.24 mg/ml

**Original Formulation:** Lyophilized powder from a protein sample in 0.1 M Glycine pH2.5, 100 mM NaCl

**Reconstitution:** Supplied as a lyophilized solid. Reconstitute in sterile dH2O.

**Stability:** ≥6 months at -80°C (as a solid). Store reconstituted protein in aliquots at -80°C for up to 6 months. Avoid freeze/thaw cycles. *Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

**References:**

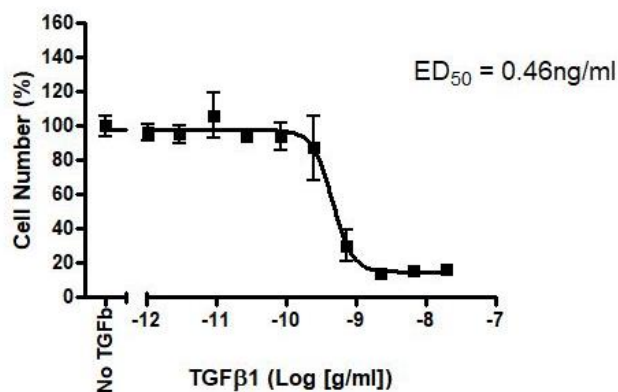
1. Derynck, R. *et al.* (1985). *Nature* **316(6030)**:701-5.
2. Bourdrel L., *et al.* (1993). *Protein Expr. Purif.* **4**:130-140.

**Description:** Human TGFβ1 (GenBank Accession #NM\_000660), a.a. 279-390(end), forms a disulfide-linked homodimer of 25 kDa (Monomer MW = 13 kDa).

**Endotoxin Level:** <1 EU/μg.

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

### Quality Assurance Specific Activity

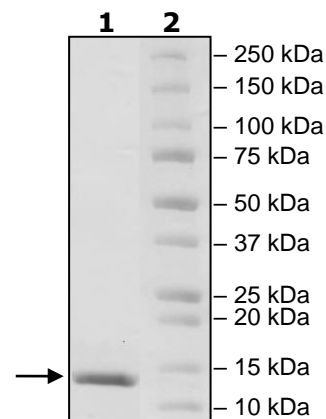


### Reducing 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2μg TGFβ1

**Lane 2:**  
Protein Marker

**MW:** 13 kDa  
**Purity:** ≥90%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)