

## Data Sheet

### TGFβ1, Latent

Human, recombinant  
**Catalog #:** 90901-3  
**Lot #:** 140813      **Size:** 1 mg

**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:** At least 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 mL-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

