

#### DESCRIPTION

**Source** *E. coli*-derived human TGF-beta 3 protein  
Accession # P10600.1

**Predicted Molecular Mass** 12.7 kDa (monomer), 25.4 kDa (dimer)

#### SPECIFICATIONS

**SDS-PAGE** Dimeric TGF- β3 protein only

**Activity** No significant difference between EC<sub>50</sub> of reference and test lots

**Endotoxin Level** <0.10 EU per 1 μg of the protein by the LAL method.

**Mass Spectrometry** Single species with expected mass

**Mycoplasma** Negative when tested in both ribosomal RNA hybridization and luminescence assays

**Formulation** Lyophilized from acetonitrile/TFA See Certificate of Analysis for details.

#### PREPARATION AND STORAGE

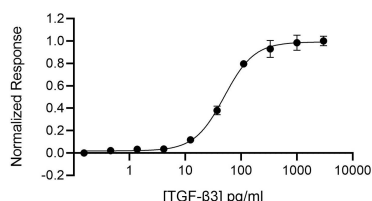
**Reconstitution** Resuspend in 10 mM HCl at >100 μg/mL, prepare single use aliquots, add carrier protein if desired.

**Shipping** The product is shipped lyophilized at ambient temperature, on ice blocks or dry ice. Shipping at ambient temperature does not affect the bioactivity or stability of the protein. Upon receipt, store immediately at the conditions stated below.

**Stability & Storage** BulkLotPrefix assignment required for Storage Info

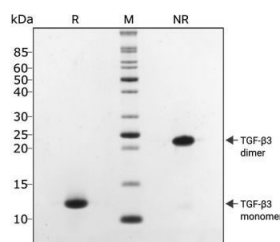
#### DATA

##### Bioactivity



**Recombinant Human TGF-beta 3, Animal-Free Protein Bioactivity** TGF-β3 activity is determined using a TGF-β3-responsive firefly luciferase reporter in HEK293T cells. Cells are treated in triplicate with a serial dilution of TGF-β3 for 6 hours. Firefly luciferase activity is measured and normalized to the control Renilla luciferase activity. EC<sub>50</sub> = 50 pg/ml (1.97 pM).

##### SDS-PAGE



**Recombinant Human TGF-beta 3, Animal-Free Protein SDS-PAGE** TGF β3 migrates as a single band at 25 kDa in non-reducing (NR) conditions and 13 kDa upon reduction (R). No contaminating protein bands are visible. Purified recombinant protein (3 μg) was resolved using 15% w/v SDS-PAGE in reduced (+β-mercaptoethanol, R) and non-reduced (NR) conditions and stained with Coomassie Brilliant Blue R250.

## BACKGROUND

TGF-β3 (transforming growth factor-beta 3) is a member of a TGF-β superfamily subgroup that is defined by their structural and functional similarities (1-5). TGF-β3 and its closely related proteins, TGF-β1 and β2, act as cellular switches to regulate immune function, cell proliferation, and epithelial-mesenchymal transition (4, 6, 7). The non-redundant biological effects of TGF-β3 include involvement in palatogenesis, chondrogenesis, and pulmonary development (1, 2, 7-9). Human TGF-β3 cDNA encodes a 412 amino acid (aa) precursor that contains a 20 aa signal peptide and a 392 aa proprotein. The proprotein is processed by a furin-like convertase to generate a 220 aa latency-associated peptide (LAP) and a 112 aa mature TGF-β3 (10, 11). Mature human TGF-β3 shows 100%, 99%, and 98% aa identity with mouse/dog/horse, rat, and pig TGF-β3, respectively. TGF-β3 is secreted as a latent complex. This latent form of TGF-β3 is activated by integrins, thrombospondin-1, plasmin, and matrix metalloproteases (12, 13). It can also be activated by extreme pH and reactive oxygen species (1-5, 12). TGF-β3 binds with high affinity to TGF-β RII, a type II serine/threonine kinase receptor. This receptor then phosphorylates and activates type I serine/threonine kinase receptors, TGF-β RI or ALK-1, to modulate transcription through Smad phosphorylation (14-16). The divergent biological effects exerted by individual TGF-β isoforms is dependent upon the recruitment of co-receptors (TGF-β RIII and endoglin) and the subsequent initiation of Smad-dependent or -independent signaling pathways (15, 17, 18).

## References:

1. Barrio, M.C. *et al.* (2014) Cells Tissues Organs. [Epub ahead of print; PMID 24861080].
2. Doetschman, T. *et al.* (2012) Genesis **50**:59.
3. Mittl, P.R. *et al.* (1996) Protein Sci. **5**:1261.
4. Sporn, M.B. (2006) Cytokine Growth Factor Rev. **17**:3.
5. Wahl, S.M. *et al.* (2006) Immunol. Rev. **213**:213.
6. Chang, H. *et al.* (2002) Endocr. Rev. **23**:787.
7. Dunker, N. and K. Kriegstein (2000) Eur. J. Biochem. **267**:6982.
8. Jin, J.Z. *et al.* (2014) Dev. Dyn. [Epub ahead of print; PMID 25104574].
9. Tang, Q.O. *et al.* (2009) Expert Opin. Biol Ther. **9**:689.
10. Derynck, R. *et al.* (1988) EMBO J. **7**:3737.
11. Miyazono, K. *et al.* (1988) J. Biol. Chem. **263**:6407.
12. Munger, J.S. *et al.* (1997) Kidney Int **51**:1376.
13. Wipff, P.J. and B. Hinz (2008) Eur J Cell Biol **87**:601.
14. Cui, X.M. and C.F. Shuler (2000) Int. J. Dev. Biol. **44**:397.
15. de Caestecker, M. (2004) Cytokine Growth Factor Rev. **15**:1.
16. Nakajima, A. *et al.* (2007) Dev. Dyn. **236**:791.
17. Iwata, J. *et al.* (2012) J. Clin. Invest. **122**:873.
18. Gatza, C.E. *et al.* (2010) Cell. Signal. **22**:1163.

## PRODUCT SPECIFIC NOTICES

The above product was manufactured, tested and released by R&D System's contract manufacturer, Qkine Ltd, at 1 Murdoch House, Cambridge, UK, CB5 8HW. The product is for research use only and not for the diagnostic or therapeutic use.