

Data Sheet

TGFβ1, Avi-Tag, Biotin-labeled

Human, recombinant, C-terminal Avi-tag,

Biotin-labeled

Catalog #: 100843

Lot#: 200807

Conc: 0.47 mg/ml

Formulated in: 0.1 M glycine, 0.1 M NaCl,
pH 2.5

Stability: At least 6 months at -80°C. *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.

Description: Human transforming growth factor beta 1, (TGFβ1), also known as TGFbeta1 and TGFB1, GenBank Accession #NM_000660, a.a. 279-390(end), with a C-terminal Avi-tag and expressed in a HEK293 expression system. MW=15 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%

Quality Assurance

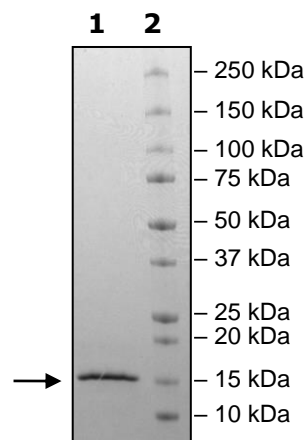
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2µg TGFbeta 1

Lane 2:
Protein Marker

MW: 15 kDa+Glycans

Purity: >90%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com