

Data Sheet

TGF β R1, Fc Fusion (IgG1) **Avi-Tag, Biotin Labeled**

Human, Recombinant, Fc-Avi-Tag,
Biotin Labeled

Catalog #: 100645

Lot #: 200409

Conc.: 0.63 mg/ml

Formulated in: 8 mM Phosphate, pH 7.4,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: At least 6 months at -80°C . Avoid
freeze/thaw cycles.

References:

1. Loeys, B.L., *et al.*, *Nature Genetics*.
2005; **37**: 275-281.
2. Singh, K.K., *et al.*, *Human Mutation*.
2006; **27(8)**: 770-777.

Description: Human Transforming growth factor beta receptor I (TGF β R1), also known as activin A receptor type II-like kinase, AAT5 and LDS1, GenBank Accession No. NM_004612.4, a.a 34-126, fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag, expressed in a HEK293 cell expression system and enzymatically biotinylated using Avi-tagTM technology. Biotinylation is confirmed to be $\geq 90\%$. MW=39 kDa. This protein runs at a higher molecular weight due to glycosylation.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

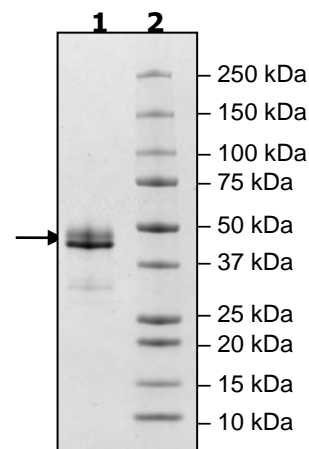
Lane 1:

4 μ g TGF β R1

Lane 2:
Protein Marker

MW: 39 kDa + glycans

Purity: $\geq 90\%$



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