

## Data Sheet

### TGF $\beta$ R2, Fc Fusion, Avi-tag, Biotin-labeled

Human, Recombinant protein,  
C-terminal Fc-Avi-Tag, Biotin-labeled  
**Catalog #:** 100629  
**Lot #:** 200227                      **Conc.:** 0.51 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  $-80^{\circ}\text{C}$ .  
Avoid freeze/thaw cycles.

#### References:

1. Lee J, *et al.*, *PLoS ONE*. 2013  
**8(2):**e57074.
2. Bhowmick, N.A, *et al.*, *Science*.  
2004; **303(5659):**848-851

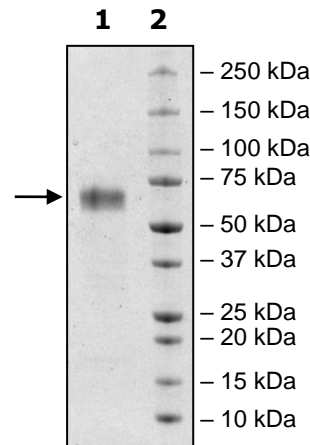
**Description:** Human transforming growth factor beta receptor 2 (TGF $\beta$ R2), also known as TGFBR2, TGFR-2, and TGF-Beta Receptor Type 2 GenBank Accession No. NM\_003242 (var2), a.a. 23-166 fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag, expressed in a HEK293 cell expression system and enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ . MW = 45 kDa. This protein runs at a higher molecular weight due to glycosylation.

## Quality Assurance

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

**Lane 1:**  
4 $\mu\text{g}$  TGF $\beta$ R2  
**Lane 2:**  
Protein Marker

**MW:** 45 kDa  
**Purity:**  $\geq 90\%$



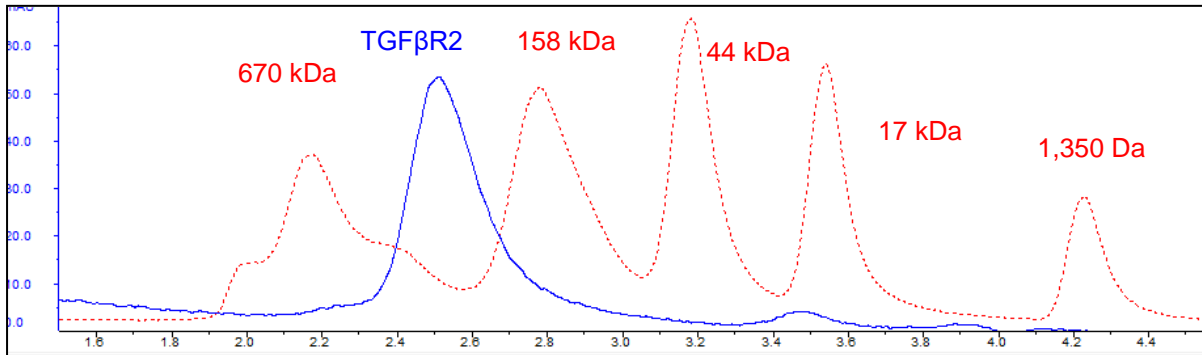
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Gel Filtration Curve



<10% Aggregation

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