

## Data Sheet

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

Mouse, Recombinant protein,  
C-terminal Fc-Avi-Tag, Biotin-labeled

**Catalog #:** 100677

**Lot #:** 200507

**Conc.:** 1.04 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. Fang, B. F., *et al.*, *Clinical & Experimental Metastasis*. 2011; **28**: 351-366
2. Sueyoshi, T. *et al.*, *Matrix Biology*. 2012; **31(6)**: 352-359.

**Description:** Mouse 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\_009371, a.a. 24-184 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 = 47 kDa. This protein runs at a higher molecular weight due to glycosylation.

---

## **Quality Assurance**

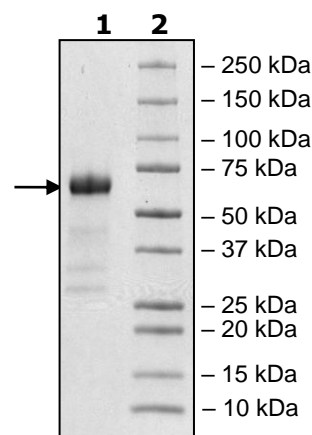
### 4-20% SDS-PAGE Coomassie staining

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

**Lane 2:**  
Protein Marker

**MW:**  
47 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](mailto:info@bpsbioscience.com)

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