

Data Sheet

TGF β R2, Fc Fusion, Avi-tag

(Mouse)

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

Catalog #: 100676

Lot #: 200408

Conc.: 1.7 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. 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 β 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, MW = 47 kDa. This protein runs at a higher molecular weight due to glycosylation.

Quality Assurance

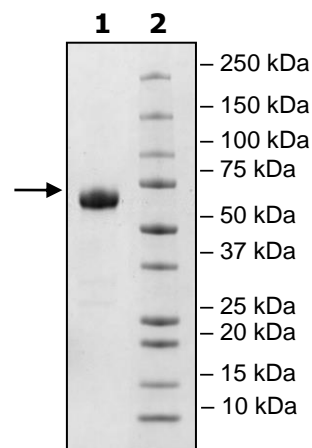
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μ g TGFbR2

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

Please visit our website at: www.bpsbioscience.com