

Data Sheet

CSF1R, Fc-Fusion, Avi-Tag (Human)

Human, Recombinant, C-terminal Fc-Avi-Tag
Catalog #: 72550
Lot #: 171010 **Conc.:** 1.52 mg/ml

Description: Human Colony stimulating factor 1 receptor (CSF1R), also known as FMS and CD115, GenBank Accession No. NM_005211, a.a. 20-512, with C-terminal Avi-Tag fused to the Fc portion of Human IgG1 and expressed in a HEK293 cell expression system. MW=83 kDa.

Formulated in: 8 mM Phosphate, pH 7.4, 400 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Walter, M., et al. (2007). *Mol. Biol.* **367** (3), 839-847
2. Li, J., et al. (2006). *Genesis* **44** (7), 328-335.

Quality Assurance

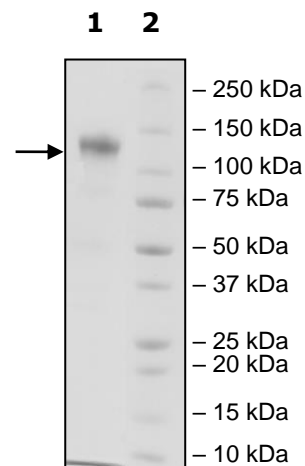
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg CD115

Lane 2:
Protein Marker

MW: 83kDa + 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 **1.858.829.3082** Fax **1.858.481.8694**

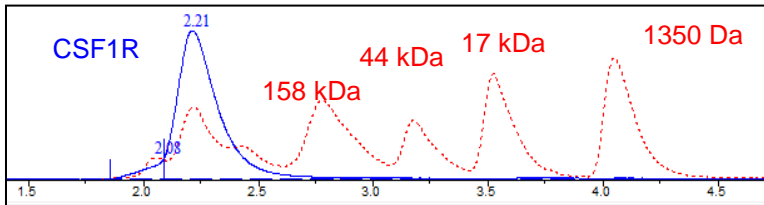
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: 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 **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com