

Data Sheet

CSF1R, Fc-Fusion, Avi-Tag, Biotin-Labeled

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

Catalog #: 79343

Lot #: 180109-G

Conc.: 0.33 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. 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.

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. This
protein is enzymatically biotinylated
using AviTag™ technology. MW = 83
kDa.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:

5µg CD115

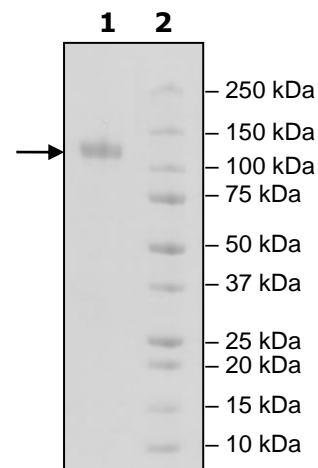
Lane 2:

Protein Marker

MW:

83kDa + glycans

Purity: ≥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