

Data Sheet

CSF1R, untagged (FMS, CD115)

Human, recombinant, N-terminal GST-tag
Catalog #: 79110

Lot #: 150616-T Conc.: 2.34 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
180 mM NaCl, 0.8 mM EDTA, 0.8 mM DTT,
0.4% Tween-20, and 25% 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. BC047521, a.a. 538-end with no tag, MW=50 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

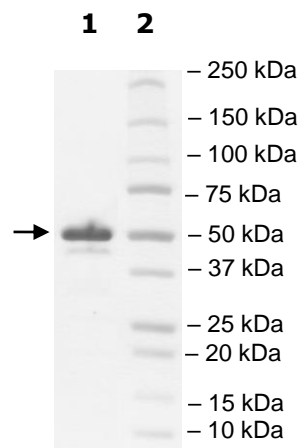
Quality Assurance

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

Lane 1:
FMS(538-e)
(2.4 ug loaded)
Lane 2:
Protein Marker

Purity: $\geq 93\%$

MW:
50 kD



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