

PRODUCT INFORMATION



CSF-1R/CD115 (human, recombinant)

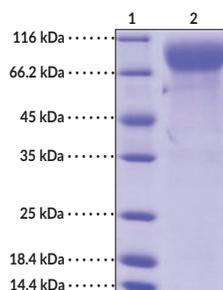
Item No. 44181

Overview and Properties

Synonyms: c-Fms, FIM2, FMS, HDLS, Macrophage Colony-stimulating Factor 1 Receptor, McDonough Feline Sarcoma Viral Oncogene Homolog, MCSF-1 Receptor
Source: Active recombinant human CSF-1R expressed in HEK293 cells
Amino Acids: 20-512
Uniprot No.: P07333
Molecular Weight: 54.5 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥90% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay
Protein Concentration: *batch specific* mg/ml

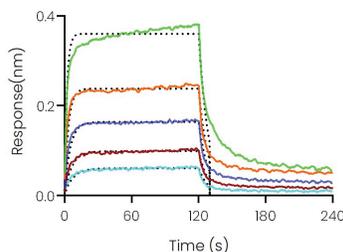
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images

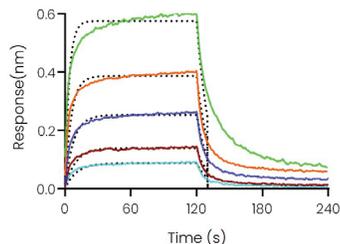


Lane 1: MW Markers
Lane 2: CSF-1R/CD115

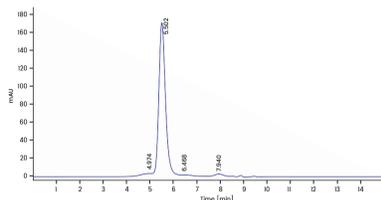
SDS-PAGE Analysis of CSF-1R/CD115. This protein has a calculated molecular weight of 54.5 kDa.



Dimerization ability of CSF-1R/CD115 in a bio-layer interferometry (BLI) assay. CSF-1R/CD115 was immobilized on a SA Biosensor via amine coupling and biotinylation. Using BLI, the affinity constant (Kd) of CSF-1R/CD115 bound to CSF-1R/CD115 was 0.323 μM.



Dimerization ability of CSF-1R/CD115 in a bio-layer interferometry (BLI) assay. CSF-1R/CD115 was immobilized on a SA Biosensor via amine coupling and biotinylation. Using BLI, the affinity constant (Kd) of CSF-1R/CD115 bound to CSF-1R/CD115 was 0.323 μM.



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/09/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Colony-stimulating factor 1 receptor (CSF-1R), also known as CD115, is a single-pass transmembrane protein encoded by *CSF1R*, previously known as the *c-fms* proto-oncogene, in humans.¹ It is composed of a glycosylated extracellular domain containing ligand-binding and dimerization regions, a transmembrane domain, and an intracellular tyrosine kinase domain. CSF-1R is expressed in monocytes, macrophages, osteoclasts, and microglia, among other cell types.² Upon binding by CSF-1R or IL-34, CSF-1R dimerizes and is autophosphorylated leading to ERK1/2, AMPK, Src/FAK, or Akt signaling.^{3,4} In this way, CSF-1R is involved in a wide variety of processes, including osteoclast cytoskeletal remodeling, macrophage motility, proliferation, and survival, and microglial proliferation.²⁻⁴ Increased intratumoral expression of *CSF1R* is associated with multiple cancer types and is predictive of poor prognosis in hematological cancers.¹ Cayman's CSF-1R/CD115 (human, recombinant) protein can be used for binding assays. This protein consists of 493 amino acids and has a calculated molecular weight of 54.5 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 100 kDa due to glycosylation.

References

1. Yeung, Y.-G. and Stanley, E.R. Proteomic approaches to the analysis of early events in colony-stimulating factor-1 signal transduction. *Mol. Cell Proteomics* **2(11)**, 1143-1155 (2003).
2. Mun, S.H., Park, P.S.U., and Park-Min, K.-H. The M-CSF receptor in osteoclasts and beyond. *Exp. Mol. Med.* **52(8)**, 1239-1254 (2020).
3. Pixley, F.J. and Stanley, E.R. CSF-1 regulation of the wandering macrophage: complexity in action. *Trends Cell Biol.* **14(11)**, 628-638 (2004).
4. Freuchet, A., Salama, A., Remy, S., et al. IL-34 and CSF-1, deciphering similarities and differences at steady state and in diseases. *J. Leukoc. Biol.* **110(4)**, 771-796 (2021).
5. Mo, H., Hao, Y., Lv, Y., et al. Overexpression of macrophage-colony stimulating factor-1 receptor as a prognostic factor for survival in cancer: A systematic review and meta-analysis. *Medicine (Baltimore)* **100(12)**, e25218 (2021).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM