

PRODUCT INFORMATION



M-CSF- β (mouse, recombinant)

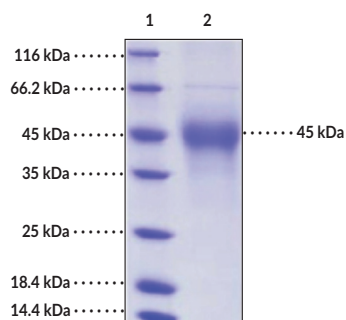
Item No. 32080

Overview and Properties

Synonym:	CSF-1
Source:	Recombinant mouse M-CSF- β expressed in HEK293 cells
Amino Acids:	33-262
Uniprot No.:	P07141-1
Molecular Weight:	26 kDa
Storage:	-80°C (as supplied)
Stability:	≥ 1 year
Purity:	$\geq 95\%$ as determined by SDS-PAGE
Supplied in:	Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing:	< 1.0 EU/ μ g, determined by the LAL endotoxin assay
Bioactivity:	Immobilized mouse M-CSF- at 10 μ g/ml (100 μ l/well) can bind mouse CSF-1R-Fch. The EC ₅₀ value of mouse CSF-1R-Fch is 0.04-0.1 μ g/ml.

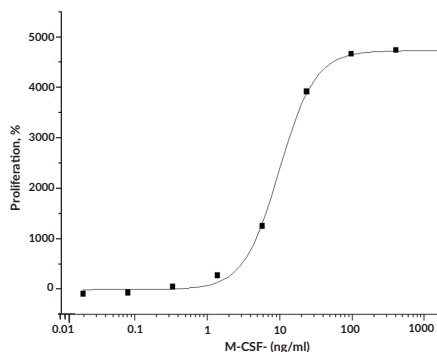
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: M-CSF- β

SDS-PAGE Analysis of M-CSF- β . This protein has a calculated molecular weight of 26 kDa. It has an apparent molecular weight of approximately 45 kDa by SDS-PAGE under reducing conditions due to glycosylation.



Cell proliferation assay using M-NFS-60 mouse myelogenous leukemia lymphoblast cells. The EC₅₀ value for this effect is typically 3-15 ng/ml.

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, 07/08/2020

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

Macrophage colony-stimulating factor (M-CSF) is a glycoprotein encoded by *Csf1* in mice that promotes the differentiation, proliferation, and function of mononuclear phagocytes, including macrophages, osteoclasts, and dendritic cells.¹ Alternative splicing of *Csf1* pre-mRNA produces one full-length long isoform (M-CSF- β), an intermediate-length isoform (M-CSF- γ), and a short-length isoform (M-CSF- α) that share sequence homology in the 150-amino acid receptor binding region which is required for the biological activity of M-CSF.^{2,3} M-CSF exists as a disulfide-linked homodimer where each monomer contains four α -helices, an anti-parallel β -sheet, and numerous glycosylation sites.¹ M-CSF is constitutively produced by many cell types, including stromal cells, osteoclasts, fibroblasts, and macrophages, and is localized to the cell surface where it can be proteolytically cleaved to yield a secreted form.^{4,5} Binding of M-CSF to the M-CSF receptor, which is expressed by monocytes, macrophages, osteoclasts, and dendritic cells, promotes cell differentiation, proliferation, and survival of mononuclear phagocytes and regulates bone resorption by osteoclasts.^{4,8} Mice homozygous for *Csf1^{op}*, an inactivating mutation, exhibit defects in fertility and neural development and develop osteopetrosis, a condition characterized by increased bone density.⁶ Neutralization of M-CSF with a monoclonal antibody decreases joint swelling and distortion in a mouse model of collagen-induced arthritis.⁷ M-CSF has been used to generate bone marrow- or peripheral blood monocyte-derived macrophages with an anti-inflammatory phenotype *in vitro*.^{8,9} Cayman's M-CSF- β (mouse, recombinant) protein can be used for binding and cell-based assay applications. This protein consists of 230 amino acids, has a calculated molecular weight of 26 kDa, and a predicted N-terminus of Lys33 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the observed molecular mass of the protein is approximately 45 kDa due to glycosylation.

References

1. Stanley, E.R., Berg, K.L., Einstein, D.B., *et al.* Biology and action of colony-stimulating factor-1. *Mol. Reprod. Dev.* **46(1)**, 4-10 (1997).
2. Cerretti, D.P., Wignall, J., Anderson, D., *et al.* Human macrophage-colony stimulating factor: Alternative RNA and protein processing from a single gene. *Mol. Immunol.* **25(8)**, 761-770 (1988).
3. Fitzgerald, K.A., Callard, R.E., O'Neill, L.A.J., *et al.* M-CSF. *The Cytokine Factsbook and Webfacts*. Fitzgerald, K.A. editor, 2nd edition., *Academic Press* (2001).
4. Chockalingam, S. and Ghosh, S.S. Macrophage colony-stimulating factor and cancer: A review. *Tumour Biol.* **35(11)**, 10635-10644 (2014).
5. Hamilton, J.A. Colony-stimulating factors in inflammation and autoimmunity. *Nat. Rev. Immunol.* **8(7)**, 533-544 (2008).
6. Dai, X.-M., Ryan, G.R., Hapel, A.J., *et al.* Targeted disruption of the mouse colony-stimulating factor 1 receptor gene results in osteopetrosis, mononuclear phagocyte deficiency, increased primitive progenitor cell frequencies, and reproductive defects. *Blood* **99(1)**, 111-120 (2002).
7. Campbell, I.K., Rich, M.J., Bischof, R.J., *et al.* The colony-stimulating factors and collagen-induced arthritis: Exacerbation of disease by M-CSF and G-CSF and requirement for endogenous M-CSF. *J. Leukoc. Biol.* **68(1)**, 144-150 (2000).
8. Stanley, E.R. and Chitu, V. CSF-1 receptor signaling in myeloid cells. *Cold Spring Harb. Mol. Case Stud.* **6(6)**, a021857 (2014).
9. Ushach, I. and Zlotnik, A. Biological role of granulocyte macrophage colony-stimulating factor (GM-CSF) and macrophage colony-stimulating factor (M-CSF) on cells of the myeloid lineage. *J. Leukoc. Biol.* **100(3)**, 481-489 (2016).

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