

# Mouse M-CSF (recombinant)



**Alias:** CSF-1 **Catalog #:** 6534

**Size:** 5 ug **Research Use Only**

**Molecular Weight:** 26.0 kDa

**Source:** Yeast. Recombinant Mouse M-CSF was produced in yeast and, therefore, does not have endotoxin. It is naturally folded and post-translationally modified.

**Formulation:** Lyophilized without carrier protein.

**Purity:** >95% as visualized by SDS-PAGE analysis.

**Purification:** Ion-exchange chromatography.

**Bioactivity:** In testing

**Entrez Gene ID:** 12977

**Number of Amino Acids:** 231

**Amino Acid Sequence:** AKEVSEHCSH MIGNGHLKVL QQLIDSQMET SCQIAFEFVD QEQLDDPVCY LKKAFFLVQD  
IIDETMRFKD NTPNANATER LQELSNLNS CFTKDYEEQN KACVRTFHET PLQLLEKIKN  
FFNETKNLLE KDWNIFTKNC NNSFAKCSSR DVVTKPDCNC LYPKATPSSD PASASPHQPP  
APSMAPLAGL AWDDSQRTEG SSSLPELPL RIEDPGSAKQ RPPRSTCQTL E

**Country of Origin:** USA

**Reconstitution:** Reconstitute with sterile phosphate-buffered saline containing at least 0.1% carrier protein.

**Stability and Storage:** Stable for up to twelve months from date of receipt at -20°C. Stable for at least 3 months when stored in working aliquots with a carrier protein at -20°C. Avoid repeated freeze/thaw cycles.

**Applications:** The M-CSF protein can be used in cell culture, as a M-CSF ELISA Standard, and as a Western Blot Control.

**Background:** Macrophage colony-stimulating factor (M-CSF), also known as CSF-1, is a hematopoietic growth factor that is involved in the proliferation, differentiation, and survival of monocytes, macrophages, and bone marrow progenitor cells. M-CSF enhances expression of differentiation antigens and stimulates chemotactic, phagocytic and the killing activities of monocytes. M-CSF also stimulates production of several cytokines, including GM-CSF, G-CSF and IL-6 by priming monocytes. It also stimulates production and secretion of IL-8 and reactive nitrogen intermediates. In addition to the stimulation of hematopoiesis, M-CSF also stimulates differentiation and proliferation of osteoclast progenitor cells and cytotrophoblasts.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Mouse M-CSF updated on: 1/29/2014