

Mouse M-CSF ELISA Kit ab155457

[2 References](#) [1 Image](#)

Overview

Product name Mouse M-CSF ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 1 pg/ml

Range 1.43 pg/ml - 350 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	98.81	93% - 106%
Plasma	103	83% - 125%
Cell culture media	133.9	127% - 138%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's Mouse M-CSF ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse M-CSF in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse M-CSF coated on a 96-well plate. Standards

and samples are pipetted into the wells and M-CSF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse M-CSF antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of M-CSF bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

Storage instructions Store at -20°C. Please refer to protocols.

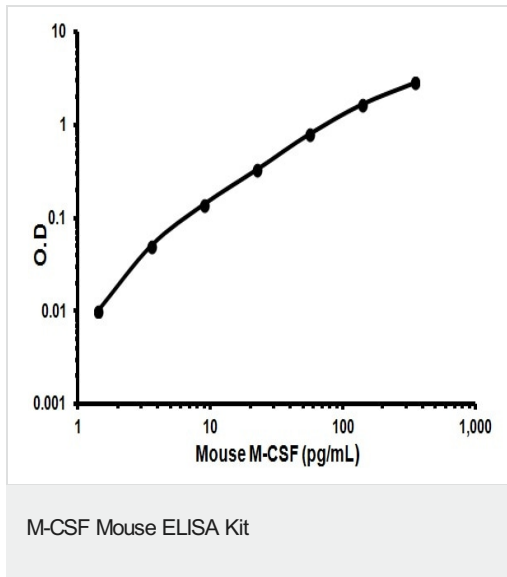
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-mouse M-CSF	2 vials
M-CSF Microplate (12 x 8 wells)	1 unit
Recombinant Mouse M-CSF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function Granulocyte/macrophage colony-stimulating factors are cytokines that act in hematopoiesis by controlling the production, differentiation, and function of 2 related white cell populations of the blood, the granulocytes and the monocytes-macrophages. CSF-1 induces cells of the monocyte/macrophage lineage. It plays a role in immunological defenses, bone metabolism, lipoproteins clearance, fertility and pregnancy.

Post-translational modifications Glycosylation and proteolytic cleavage yield different soluble forms. A high molecular weight soluble form is a proteoglycan containing chondroitin sulfate. Isoform 1 is N- and O-glycosylated. Isoform 3 is N-glycosylated.

Cellular localization Cell membrane and Secreted > extracellular space.

Images



Representative standard curve using ab155457

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors