

Mouse CSF-1-R ELISA Kit ab213858

1 Image

Overview

Product name Mouse CSF-1-R ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	106pg/ml	8.16	= 7.7%
2	16	398pg/ml	24.27	= 6.1%
3	16	1799pg/ml	131.32	= 7.3%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	10.23pg/ml	10.23	= 9.3%
2	24	32.87pg/ml	32.87	= 7.7%
3	24	137.74pg/ml	137.74	= 7.8%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 62.5 pg/ml - 4000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview The Mouse CSF-1-R ELISA Kit (ab213858) is designed for the quantitative measurement of mouse CSF-1-R in cell culture supernatants, cell lysates, serum or plasma (heparin, EDTA).

It is based on standard sandwich enzyme-linked immunosorbent assay technology. A monoclonal antibody from rat specific for CSF-1-R has been pre-coated onto 96-well plates. Standards

(Expression system for standard: NSO; Immunogen sequence: A20-S511) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for CSF-1-R is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the mouse CSF-1-R amount of sample captured in plate.

Notes CSF-1-R (Colony stimulating factor 1 receptor), also known as M-CSF-R and CD115, is a cell-surface protein encoded, in humans, by the CSF1R gene. The gene is located on long arm of chromosome 5 (5q32) on the Crick (minus) strand. The encoded protein is a tyrosine kinase transmembrane receptor and member of the CSF1/PDGF receptor family of tyrosine-protein kinases. The encoded protein is a single pass type I membrane protein and acts as the receptor for colony stimulating factor 1, a cytokine which controls the production, differentiation, and function of macrophages. Both CSF-1-R and its ligand colony stimulating factor 1 play an important role in the development of the mammary gland and may be involved in the process of mammary gland carcinogenesis.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Mouse CSF-1-R coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Mouse CSF-1-R antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Mouse CSF-1-R standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function Protein tyrosine-kinase transmembrane receptor for CSF1 and IL34.

Tissue specificity Expressed in bone marrow and in differentiated blood mononuclear cells.

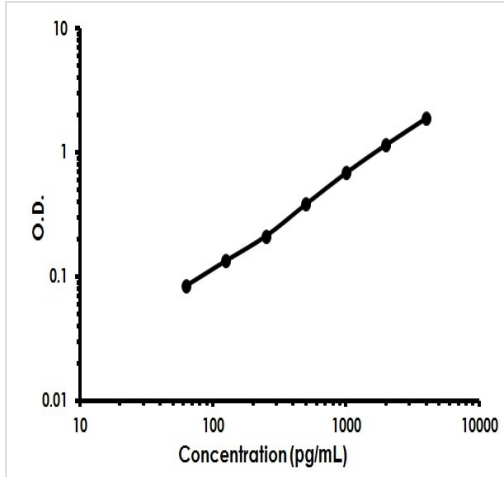
Sequence similarities Belongs to the protein kinase superfamily. Tyr protein kinase family. CSF-1/PDGF receptor subfamily.

Contains 5 Ig-like C2-type (immunoglobulin-like) domains.
Contains 1 protein kinase domain.

Cellular localization

Membrane.

Images



Mouse CSF-1-R ELISA Kit (ab213858) Standard Curve.

Mouse CSF-1-R ELISA Kit (ab213858) Standard
Curve

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