

Mouse G-CSF ELISA Kit (CSF3) ab100684

2 Images

Overview

Product name	Mouse G-CSF ELISA Kit (CSF3)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.5 pg/ml
Range	0.62 pg/ml - 150 pg/ml
Recovery	95 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	97.52	84% - 104%
Serum	95.53	83% - 104%
Plasma	94.28	82% - 102%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Mouse

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

This assay employs an antibody specific for mouse G-CSF coated on a 96- well plate. Standards and samples are pipetted into the wells and G-CSF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse G-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 G-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

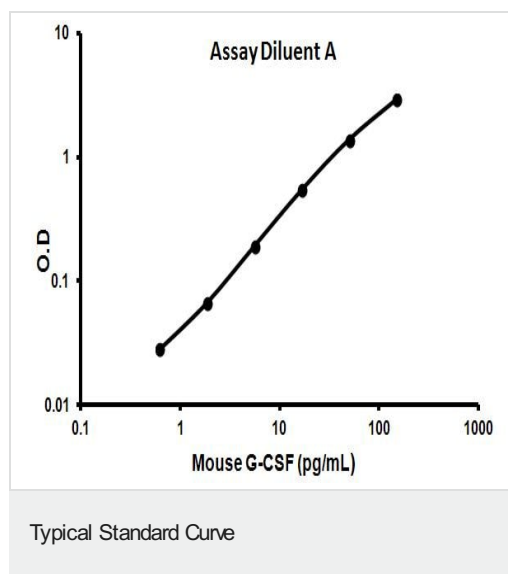
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
80x HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse G-CSF	2 vials
G-CSF Microplate (12 x 8 wells)	1 unit
Recombinant Mouse G-CSF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

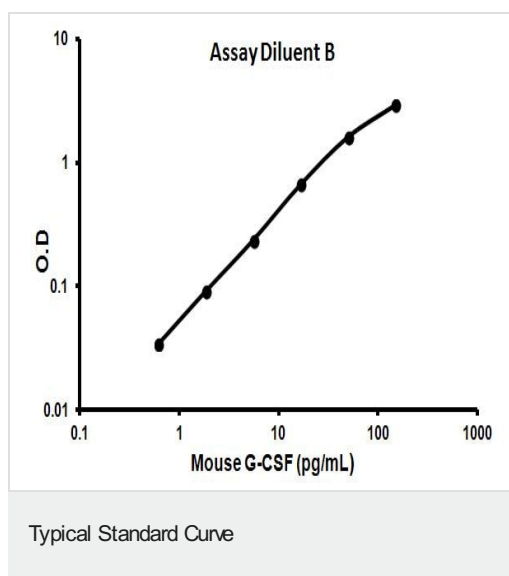
Relevance

Granulocyte Colony Stimulating Factor (G-CSF) a cytokine that controls the production, differentiation, and function of granulocytes. It is a potent stimulator of bone marrow cells especially those of neutrophil lineage. In addition, G-CSF can enhance the survival and activate the immunological functions of mature neutrophils. Recombinant human G-CSF has been developed by pharmaceutical companies, and since the late 1980's, multiple clinical trials have explored its efficacy in a variety of conditions.

Images



Representative Standard Curve using ab100684



Representative Standard Curve using ab100684

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