

# Human G-CSF ELISA Kit ab100524

[2 References](#)   [5 Images](#)

Overview

Product name	Human G-CSF ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1 pg/ml
Range	0.69 pg/ml - 500 pg/ml
Recovery	> 75 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	76.56	68% - 90%
Serum	83.45	72% - 94%
Plasma	74.76	67% - 87%

Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Human
Product overview	Abcam's G-CSF Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human G-CSF in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human 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-Human 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.

Notes	Optimization may be required with urine samples
Platform	Microplate

## Properties

### Storage instructions

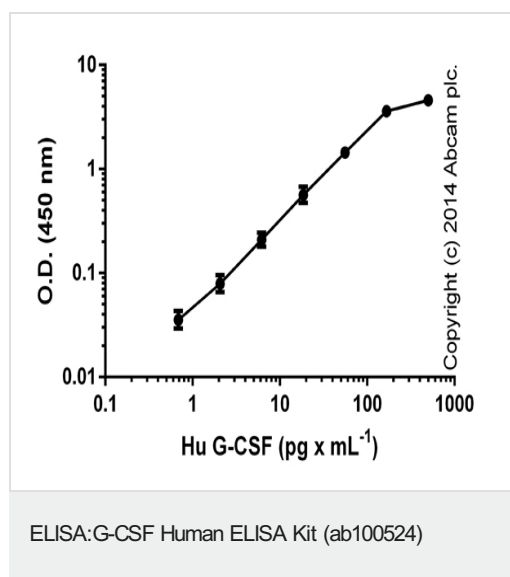
Store at -20°C. Please refer to protocols.

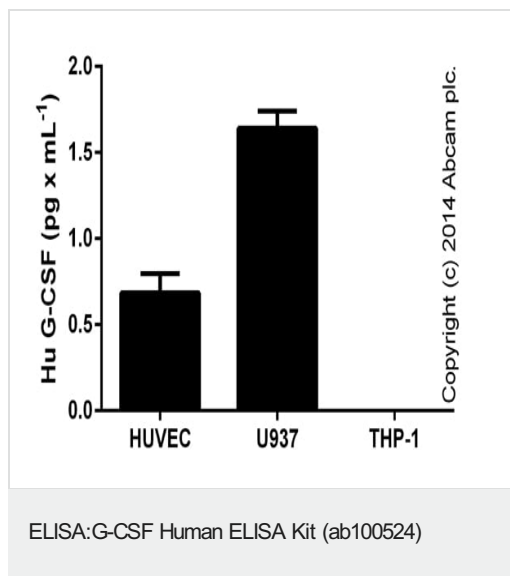
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
260X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human G-CSF	2 vials
Human G-CSF Microplate (12 x 8 wells)	1 unit
Recombinant Human G-CSF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### Relevance

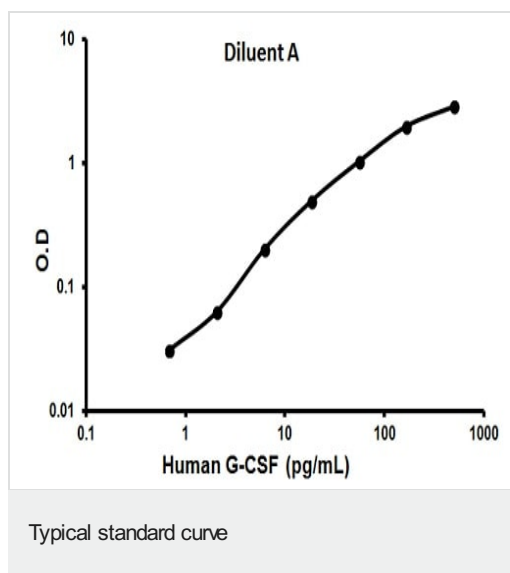
Granulocyte Colony Stimulating Factor (G-CSF) a cytokine that controls the production, differentiation, and function of granulocytes. It is a potent stimulator of bone marrow cells especially those of neutrophil lineage. In addition, G-CSF can enhance the survival and activate the immunological functions 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

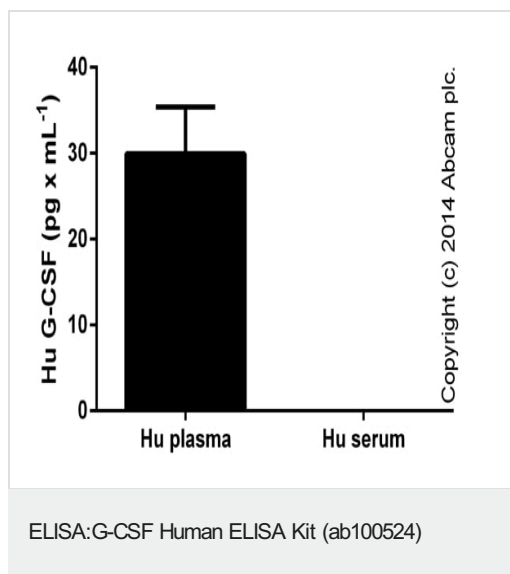




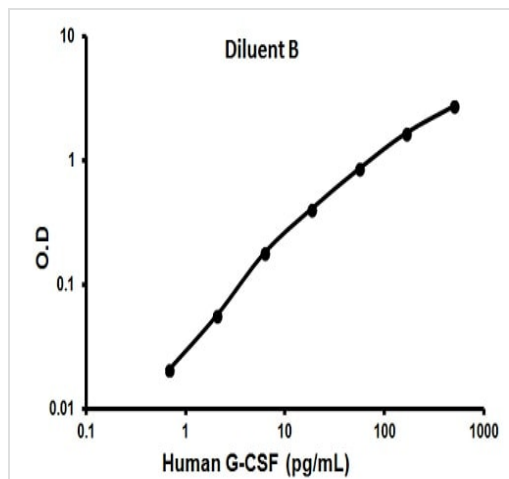
Human G-CSF measured in undiluted culture supernatants (duplicates +/- SD).



Representative standard curve using ab100524



G-CSF measured in undiluted human plasma, signal in serum was below level of detection (duplicates +/- SD).



Representative standard curve using ab100524

Typical 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