

Human G-CSF ELISA Kit ab100524

画像数 5

製品の概要

製品名	Human G-CSF ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 1 pg/ml
検出範囲	0.69 pg/ml - 500 pg/ml
添加回収試験	> 75 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	76.56	68% - 90%
Serum	83.45	72% - 94%
Plasma	74.76	67% - 87%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Abcam's G-CSF Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* 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.

Get results in 90 minutes with Human G-CSF ELISA Kit (CSF3) ([ab188390](#)) from our

SimpleStep ELISA[™] range.

特記事項	Optimization may be required with urine samples
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Microplate

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer Concentrate	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 (lyophilized)	2 vials
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

関連性 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.

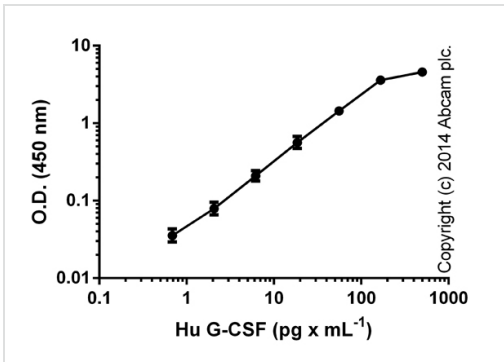
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab100524** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

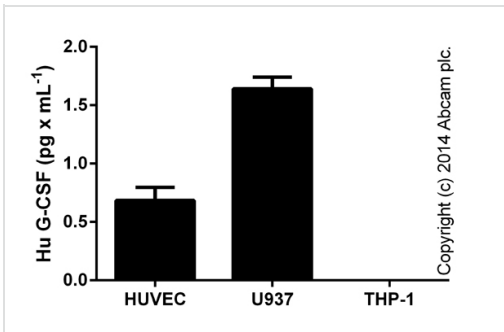
アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



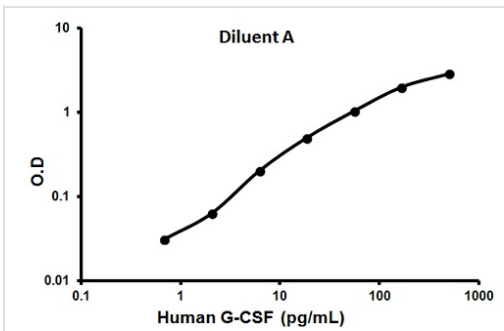
Standard curve in diluent B with background signal subtracted (duplicates; +/- SD).

ELISA:G-CSF Human ELISA Kit (ab100524)



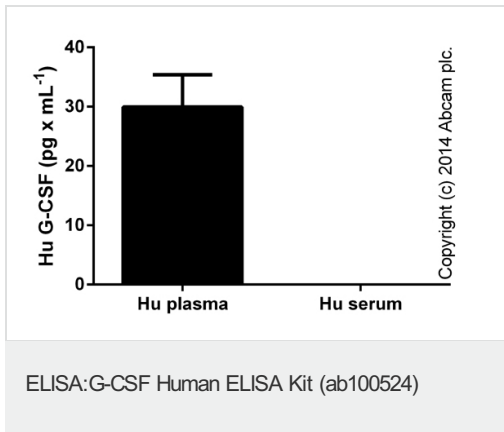
Human G-CSF measured in undiluted culture supernatants (duplicates +/- SD).

ELISA:G-CSF Human ELISA Kit (ab100524)

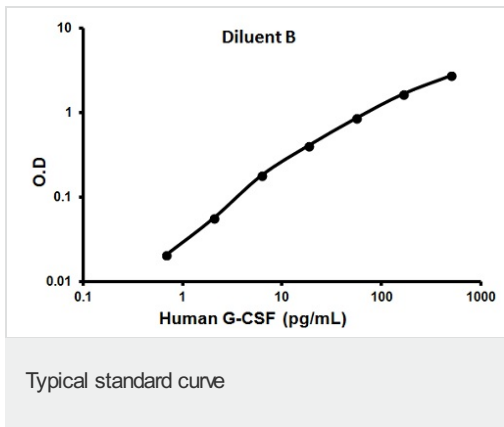


Representative standard curve using ab100524

Typical standard curve



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



Representative standard curve using ab100524

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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