

Human G-CSF R ELISA Kit ab267649

[1 Image](#)

Overview

Product name Human G-CSF R 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 0.1 ng/ml

Range 0.1 ng/ml - 25 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	86.34	70% - 99%
Serum	81.01	69% - 96%
Plasma	77.54	68% - 92%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human G-CSF R ELISA Kit is designed for the quantitative determination of G-CSF R in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human G-CSF R coated on a 96-well plate. Standards and samples are pipetted into the wells and G-CSF R present in a sample is bound to

the wells by the immobilized antibody. The wells are washed, and biotinylated anti-human G-CSF R 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 R bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit shows no cross-reactivity with the following cytokines tested: Human ACE-2, aFGF (FGF-1), Albumin, AMICA, Angiopoietin-4, BAFF, CA19-9, CD163, Clusterin, CRTAM, CXCL14, Cystatin C, Decorin, Dkk-3, DLL1, Fetuin A, FOLR1, Furin, GASP-1, GASP-2, HAI-2, IL-17 RB, IL-27, LAG-3, LDL R, Pepsinogen 1, RANK, RBP4, SOST, Syndecan-1, TACI, TFPI, Thrombospondin 1, TRAIL R2, TRANCE, Troponin I, uPA, VE-Cadherin, WISP-1.

Platform Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Anti-Human G-CSF R coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human G-CSF R Detection Antibody	2 vials
Human G-CSF R Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Receptor for granulocyte colony-stimulating factor (CSF3), essential for granulocytic maturation. Plays a crucial role in the proliferation, differentiation and survival of cells along the neutrophilic lineage. In addition it may function in some adhesion or recognition events at the cell surface.

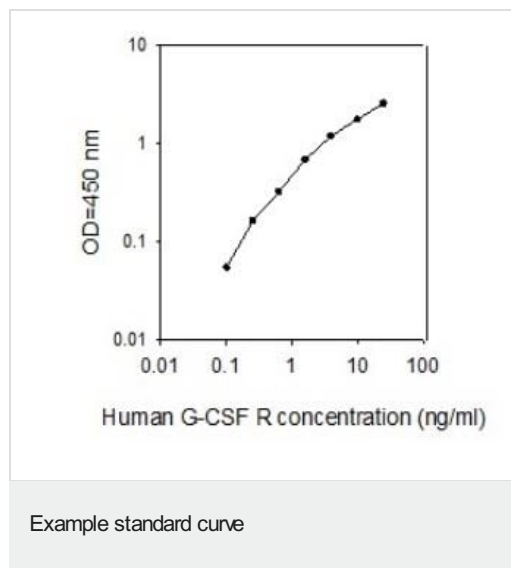
Tissue specificity One or several isoforms have been found in myelogenous leukemia cell line KG-1, leukemia U-937 cell line, in bone marrow cells, placenta, and peripheral blood granulocytes. Isoform GCSFR-2 is found only in leukemia U-937 cells. Isoform GCSFR-3 is highly expressed in placenta.

Involvement in disease Hereditary neutrophilia

Sequence similarities Belongs to the type I cytokine receptor family. Type 2 subfamily.
Contains 5 fibronectin type-III domains.
Contains 1 Ig-like C2-type (immunoglobulin-like) domain.

Domain The WSXWS motif appears to be necessary for proper protein folding and thereby efficient intracellular transport and cell-surface receptor binding.
The box 1 motif is required for JAK interaction and/or activation.

Images



Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.

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