

## Product datasheet

# Human GM-CSF ELISA Kit ab100529

1 Image

### Overview

<b>Product name</b>	Human GM-CSF ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 2 pg/ml
<b>Range</b>	2.05 pg/ml - 500 pg/ml
<b>Recovery</b>	> 75 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	101.1	89% - 114%
Serum	75.43	69% - 83%
Plasma	74.49	68% - 83%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

### Product overview

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

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

**Tested applications** **Suitable for:** Sandwich ELISA

**Platform** Microplate

## Properties

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

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

**Function** Cytokine that stimulates the growth and differentiation of hematopoietic precursor cells from various lineages, including granulocytes, macrophages, eosinophils and erythrocytes.

**Sequence similarities** Belongs to the GM-CSF family.

**Cellular localization** Secreted.

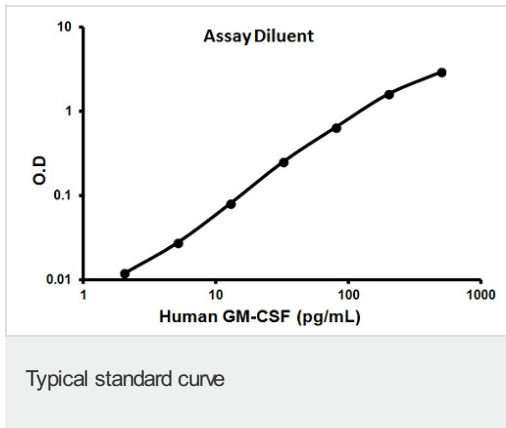
## Applications

Our [Abpromise guarantee](#) covers the use of **ab100529** in the following tested applications.

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

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



**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 <http://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