

Human GM-CSF Receptor alpha ELISA Kit (CD116)
ab46086

1 Image

Overview

Product name	Human GM-CSF Receptor alpha ELISA Kit (CD116)																																	
Detection method	Colorimetric																																	
Precision	<div>Intra-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>A</td><td>31</td><td>700.2pg/ml</td><td>20.4</td><td>2.9%</td></tr><tr><td>B</td><td>32</td><td>378pg/ml</td><td>12.8</td><td>3.4%</td></tr></table> <div>Inter-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>A</td><td>6</td><td>555pg/ml</td><td>22.2</td><td>4%</td></tr><tr><td>B</td><td>6</td><td>120.7pg/ml</td><td>2.2</td><td>1.8%</td></tr></table>				Sample	n	Mean	SD	CV%	A	31	700.2pg/ml	20.4	2.9%	B	32	378pg/ml	12.8	3.4%	Sample	n	Mean	SD	CV%	A	6	555pg/ml	22.2	4%	B	6	120.7pg/ml	2.2	1.8%
Sample	n	Mean	SD	CV%																														
A	31	700.2pg/ml	20.4	2.9%																														
B	32	378pg/ml	12.8	3.4%																														
Sample	n	Mean	SD	CV%																														
A	6	555pg/ml	22.2	4%																														
B	6	120.7pg/ml	2.2	1.8%																														
Sample type	Cell culture supernatant, Serum, Plasma																																	
Assay type	Sandwich (quantitative)																																	
Sensitivity	< 16 pg/ml																																	
Range	31.25 pg/ml - 1000 pg/ml																																	
Assay time	3h 45m																																	
Assay duration	Multiple steps standard assay																																	
Species reactivity	Reacts with: Human																																	
Product overview	<p>Abcam's GM-CSF Receptor alpha (CD116) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of GM-CSF R alpha in serum, plasma, and cell culture supernatants.</p> <p>A monoclonal antibody specific for GM-CSF R alpha has been coated onto the wells of the microtiter strips provided. Samples, including standards of known GM-CSF R alpha concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for GM-</p>																																	

CSF R alpha are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of GM-CSF R alpha present in the samples.

This kit will recognize both endogenous and recombinant Human GM-CSF R alpha.

Platform

Microplate

Properties

Storage instructions

Store at +4°C. Please refer to protocols.

Components	Identifier	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Biotinylated anti-GM-CSF R alpha	Red	2 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
Control	Silver	4 vials	2 vials
GM-CSF Receptor alpha Microplate (12 x 8 well strips)		2 units	1 unit
GM-CSF Receptor alpha Standard (Lyophilized)	Yellow	4 vials	2 vials
HRP Diluent	Red	1 x 23ml	1 x 12ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

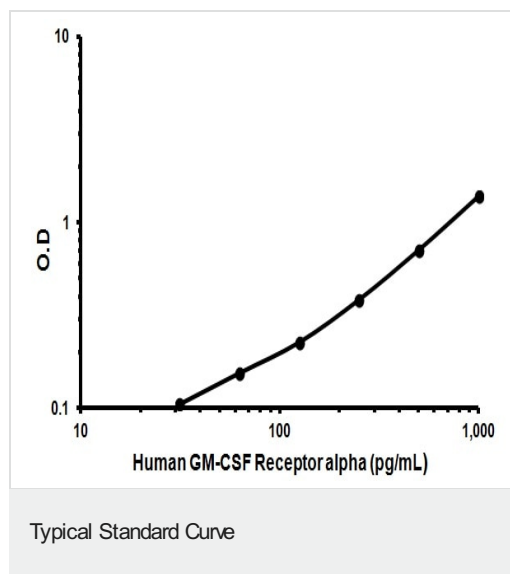
Relevance

GM-CSF Receptor alpha is the alpha subunit of the heterodimeric receptor for colony stimulating factor 2, a cytokine which controls the production, differentiation, and function of granulocytes and macrophages. The encoded protein is a member of the cytokine family of receptors. This gene is found in the pseudoautosomal region of the X and Y chromosomes. Multiple transcript variants encoding different isoforms have been found for this gene, with some of the isoforms being membrane-bound and others being soluble.

Cellular localization

Cell membrane; single-pass type I membrane protein. Isoform 3, isoform 4, isoform 6: Secreted protein

Images



Representative Standard Curve using ab46086

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