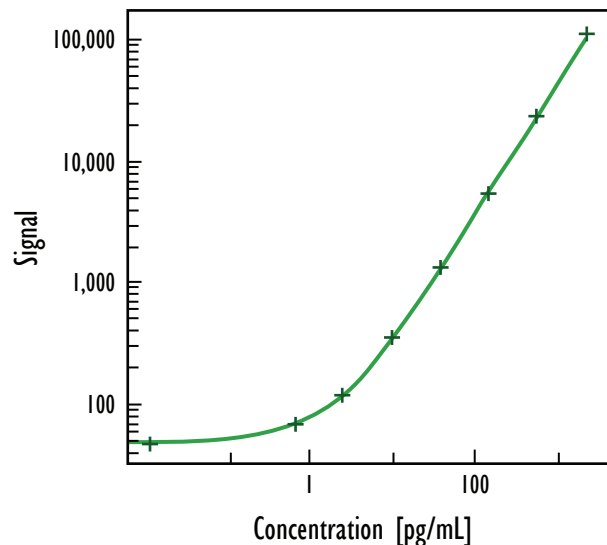


MULTI-ARRAY® Human Granulocyte Colony Stimulating Factor Ultra-Sensitive Assay

Detection of Granulocyte Colony Stimulating Factor (G-CSF) in Human Serum and Plasma Samples

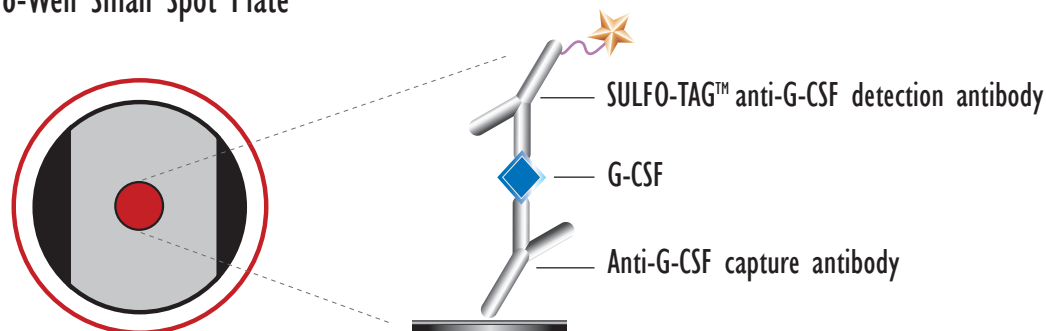
Typical Standard Curve Data



Human G-CSF	
Concentration (pg/mL)	Mean Signal
0	48
0.6	70
2.4	123
10	356
39	1,364
156	5,346
625	23,344
2,500	115,151

Standard curve data is from a representative experiment

MSD MULTI-ARRAY®
96-Well Small Spot Plate



Typical Detection Parameters

Protocol	LLOD (pg/mL)
Serum and Plasma	1.5

Definition of Detection Parameters

LLOD = Lower Limit of Detection is
2.5 stdev above the background

Kit Size	Catalog Number
1 plate	K151IPC-1
5 plates	K151IPC-2
20 plates	K151IPC-3
20 plates (Base)	K151IPC-3

MULTI-ARRAY® Human Granulocyte Colony Stimulating Factor Ultra-Sensitive Assay

Detection of Granulocyte Colony Stimulating Factor (G-CSF) in Human Serum and Plasma Samples

Spike Recovery

	Spike Level (pg/mL)	% Recovery of Spiked Calibrator
Serum	39	79%
	156	87%
EDTA Plasma	39	77%
	156	78%
Heparin Plasma	39	72%
	156	73%

- Measured analyte spiked into neat human samples
- % recovery = $\frac{\text{measured value} * 100}{\text{expected value}}$

Endogenous Levels of Human G-CSF

	Endogenous Analyte Levels, pg/mL	
Serum	Mean	15
	CV	2%
EDTA Plasma	Mean	13
	CV	11%
Heparin Plasma	Mean	11
	CV	11%

- Pooled normal human serum and plasma samples were tested for endogenous G-CSF
- Detected level was above LOQ for each sample type

Dilution Linearity

	Dilution Factor	% Recovery of Dilution Linearity
Serum	1/2	118%
	1/4	114%
	1/8	< LOQ
EDTA Plasma	1/2	125%
	1/4	130%
	1/8	< LOQ
Heparin Plasma	1/2	116%
	1/4	132%
	1/8	< LOQ

- Serum and plasma samples were diluted in HSC Assay Diluent prior to assay
- % recovery = $\frac{\text{measured value} * \text{dilution factor} * 100}{\text{predicted value}}$
- Dilutions that resulted in counts values less than the lower limit of quantitation are shown as "< LOQ"