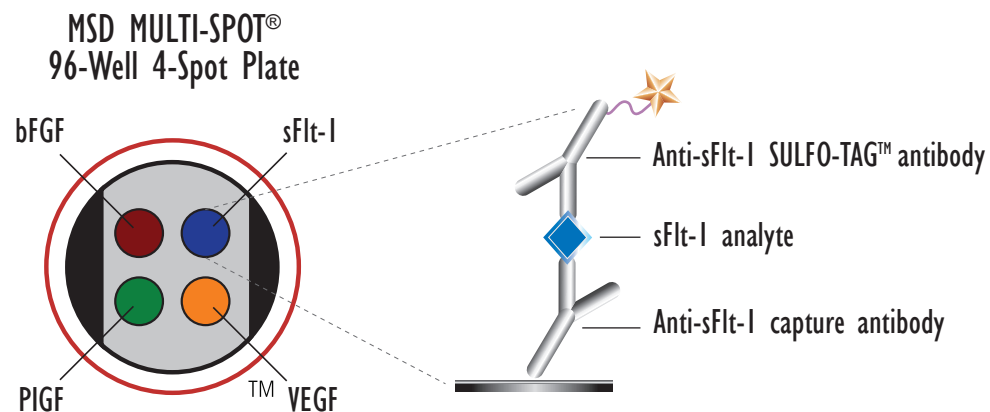
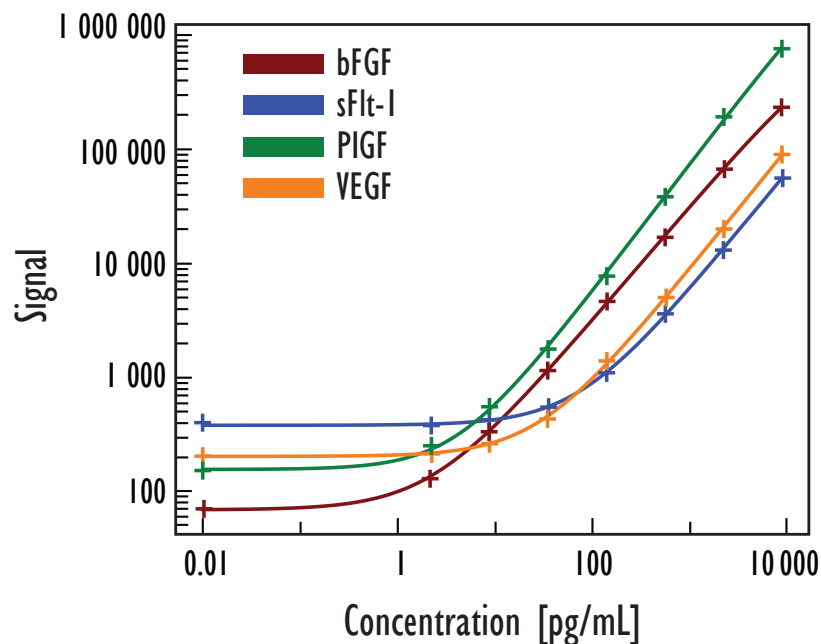


# MULTI-SPOT<sup>®</sup> Human Growth Factor Panel I Assay

Detection of bFGF, sFlt-1, PlGF, and VEGF in Human Serum and Plasma Samples



	bFGF	sFlt-1	PlGF	VEGF
LLOD (pg/mL):	2.2	8.9	0.96	6.4

LLOD (Lower Limit of Detection) is defined as 2.5x stdev above the background.

Calibrator (pg/mL)	bFGF		sFlt-1		PlGF		VEGF	
	Signal	%CV	Signal	%CV	Signal	%CV	Signal	%CV
0	71	4.2	404	4.3	150	8.3	208	4.8
2	133	3.7	387	7.6	250	4.7	218	7.1
9	345	2.6	415	3.0	558	0.3	265	6.1
35	1149	5.4	543	3.3	1784	4.0	437	4.6
141	4641	1.7	1129	2.4	7928	1.1	1422	0.8
563	17326	5.0	3550	3.4	38257	3.4	4903	2.5
2250	65545	3.1	13238	3.6	190447	3.2	20087	2.5
9000	235652	2.2	55940	3.9	765973	3.0	90165	0.4

Above is representative calibration curve data

Kit Size	Catalog Number
1 plate	K15029C-1
5 plates	K15029C-2
20 plates	K15029C-3
20 plates (Base)	K15029A-3

# MULTI-SPOT<sup>®</sup> Human Growth Factor Panel I Assay

*Detection of bFGF, sFlt-1, PlGF, and VEGF in Human Serum and Plasma Samples*

## Recovery of Spiked Calibrator

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

	Spike Level (pg/mL)	Serum	EDTA Plasma
bFGF	100	99	77
sFlt-1	1,000	105	94
PlGF	100	111	98
VEGF	100	114	95

## Recovery of Dilution Linearity

- Measured analyte levels in pooled serum diluted in assay diluent
- % recovery =  $\frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$

Dilution Factor	bFGF	sFlt-1	PlGF	VEGF
0.75	107	107	104	105
0.5	106	101	101	99
0.25	104	98	98	93

## Endogenous Analyte Levels in Samples

- Normal pooled samples

	bFGF (pg/mL)	sFlt-1 (pg/mL)	PlGF (pg/mL)	VEGF (pg/mL)
Serum	< 1	63	13	31
EDTA Plasma	7	201	16	148