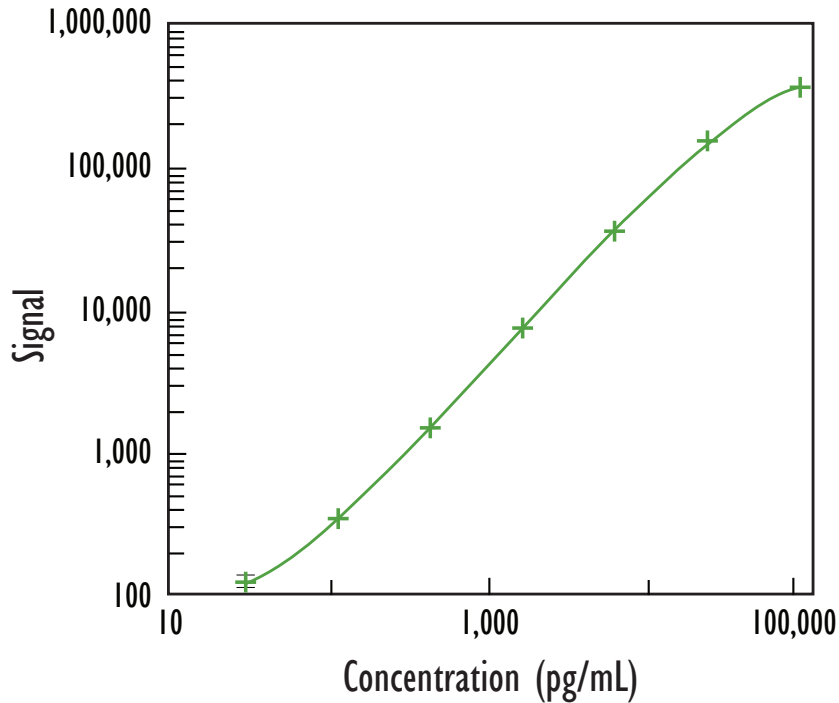


MULTI-ARRAY[®] Human IGFBP-1 Tissue Culture Assay

Detection of Human Insulin-like Growth Factor Binding Protein-1 (IGFBP-1) in Tissue Culture

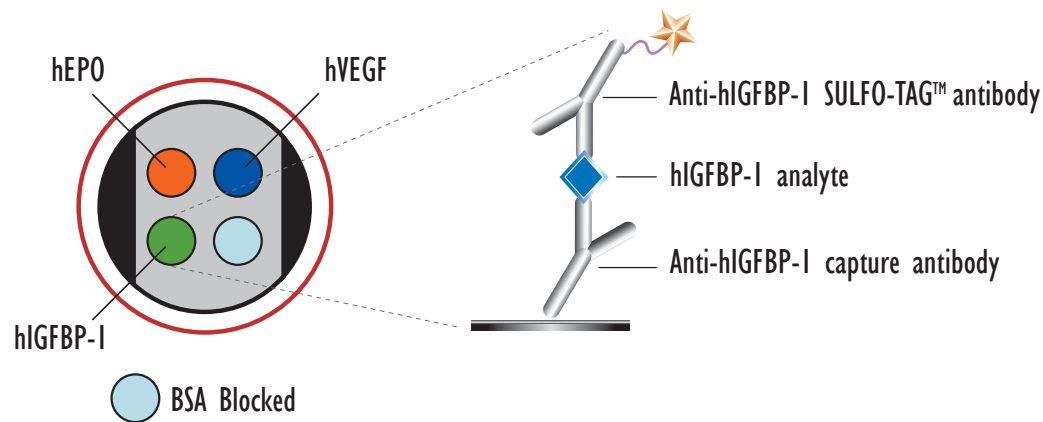


Concentration (pg/mL)	hIGFBP	
	Mean	%CV
0	102	16
24	205	10
98	512	3
391	1,872	3
1,563	7,901	3
6,250	32,446	1
25,000	118,220	2
100,000	256,565	1

Above is representative calibration curve data.

	hIGFBP
Average LLOD	15 pg/mL

LLOD is defined as 2.5x stdev above the background.



Kit Size	Catalog Number
1 plate	K151BRB-1
5 plates	K151BRB-2
20 plates	K151BRB-3
20 plates (Base)	K151BRA-3