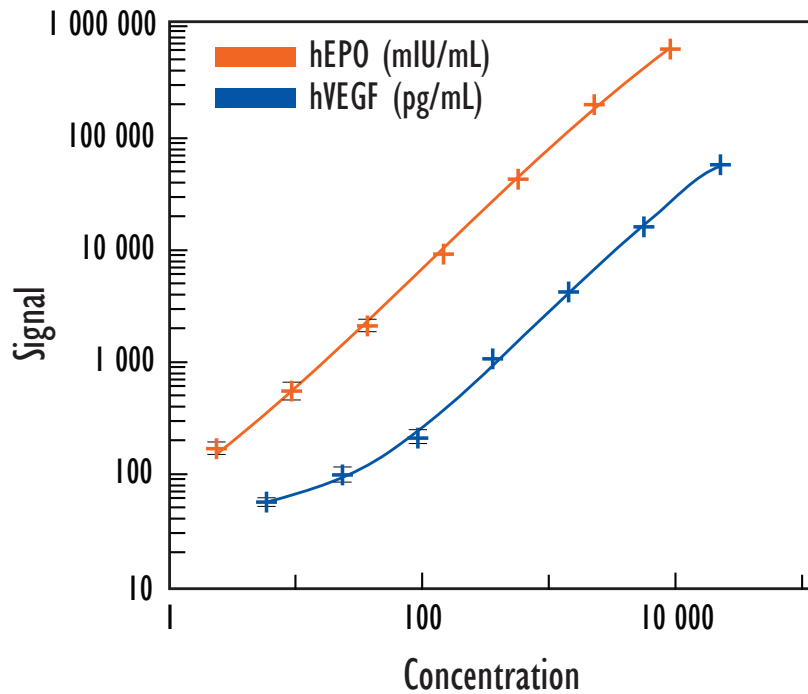


MULTI-SPOT[®] Human Hypoxia Tissue Culture Assay

Detection of Human Erythropoietin (EPO) and Vascular Endothelial Growth Factor in Tissue Culture

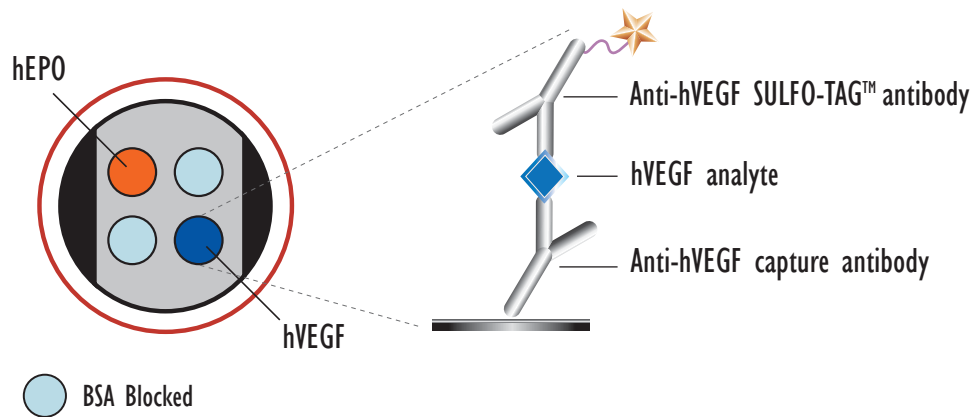


Concentration (mIU/mL)	hEPO		Concentration (pg/mL)	hVEGF	
	Mean	%CV		Mean	%CV
0	48	47	0	48	30
2	171	13	6	56	10
10	554	19	24	100	15
39	2 010	7	98	217	15
156	9 190	3	391	1 066	3
625	42 080	4	1 563	4 212	4
2 500	196 083	2	6 250	14 922	9
10 000	611 690	2	25 000	56 854	3

Above is representative calibration curve data.

	hEPO	hVEGF
Average LLOD	1 mIU/mL	9 pg/mL

LLOD is defined as 2.5x stdev above the background.



Kit Size	Catalog Number
1 plate	K15122B-1
5 plates	K15122B-2
20 plates	K15122B-3
20 plates (Base)	K15122A-3