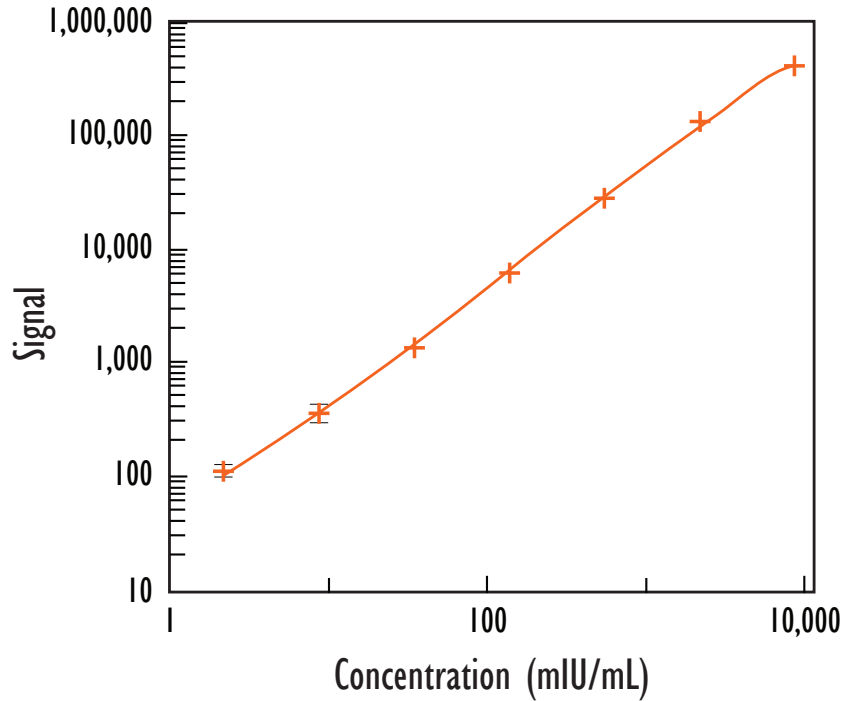


# MULTI-ARRAY<sup>®</sup> Human EPO Tissue Culture Assay

## Detection of Human Erythropoietin (EPO) in Tissue Culture

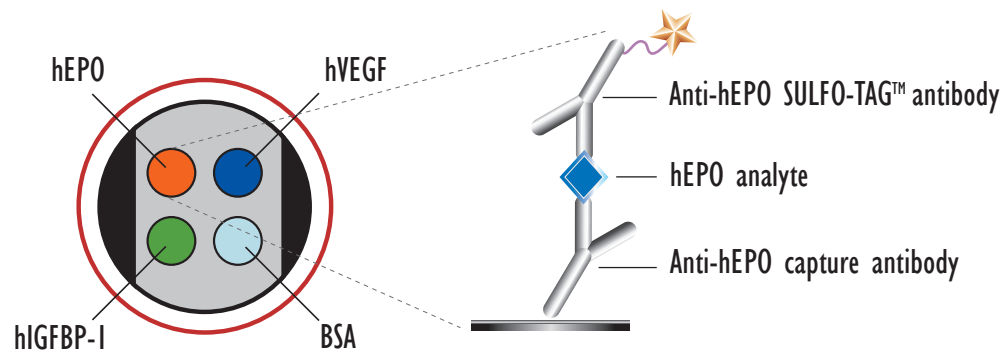


Concentration (mIU/mL)	hEPO	
	Mean	%CV
0	48	47
2	171	13
10	554	19
39	2,010	7
156	9,190	3
625	42,080	4
2,500	196,083	2
10,000	611,690	2

Above is representative calibration curve data.

	hEPO
Average LLOD	1 mIU/mL

LLOD is defined as 2.5x stdev above the background.



Kit Size	Catalog Number
1 plate	K151BQB-1
5 plates	K151BQB-2
20 plates	K151BQB-3
20 plates (Base)	K151BQA-3