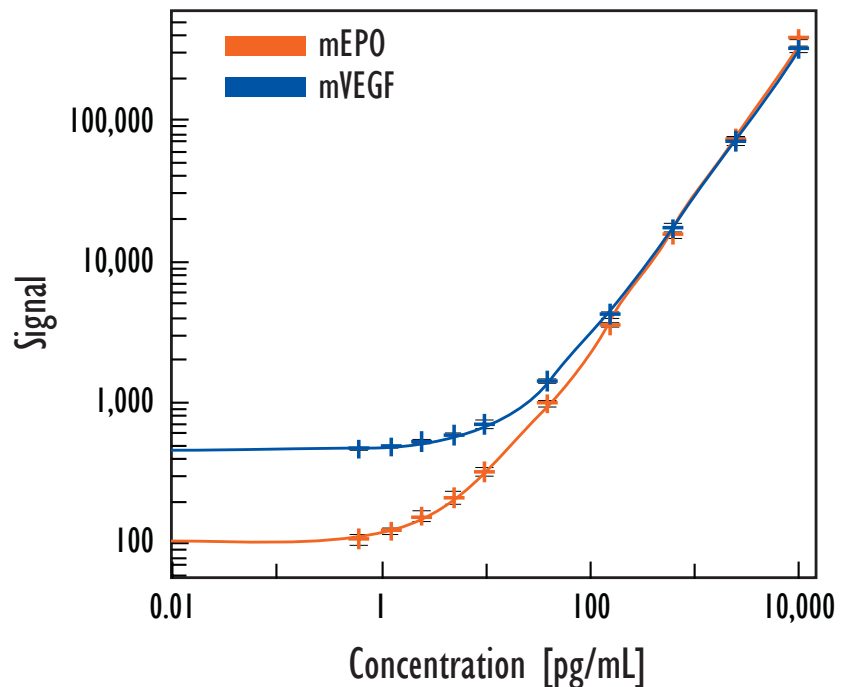


# MULTI-SPOT<sup>®</sup> Mouse/Rat Hypoxia Tissue Culture Kit

*Assay Performance for Mouse Samples Using the Tissue Culture Protocol*



Concentration (pg/mL)	mEPO		mVEGF	
	Mean	%CV	Mean	%CV
0	100	12.4	449	3.5
0.61	107	7.9	467	2.8
1.2	123	4.9	485	2.7
2.4	156	7.8	529	2.4
4.9	213	9.2	593	1.7
10	325	7.4	697	6.2
39	979	5.3	1,426	3.2
156	3,545	4.7	4,213	5.4
625	15,305	5.4	17,139	8.4
2,500	74,237	0.9	70,988	7.4
10,000	380,026	0.8	319,323	6.0

Standard curve data is from a representative experiment.

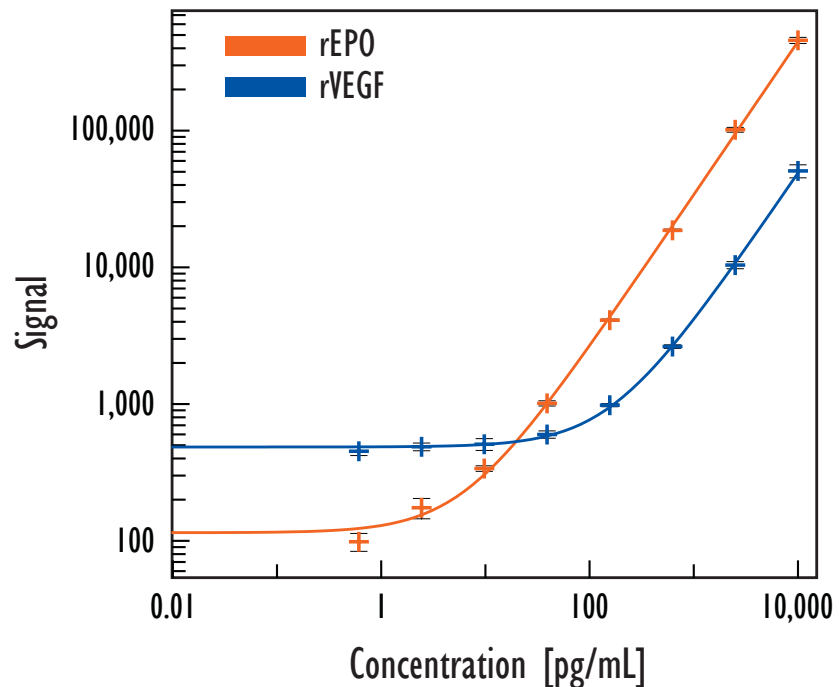
Kit Size	SI2400	SI6000
1 plate	K15123B-1	K11123B-1
5 plates	K15123B-2	K11123B-2
20 plates	K15123B-3	K11123B-3
20 plates (Base)	K15123A-3	K11123A-3

Detection Limit (pg/mL)	mEPO	mVEGF
Average LLOD	1.4	2.1

LLOD is defined as 2.5x stdev above the background.

# MULTI-SPOT<sup>®</sup> Mouse/Rat Hypoxia Tissue Culture Kit

## Assay Performance for Rat Samples Using the Tissue Culture Protocol



Concentration (pg/mL)	rEPO		rVEGF	
	Mean	%CV	Mean	%CV
0	152	14	514	7
0.6	99	15	452	7
2	175	17	487	6
10	338	5	509	10
39	1,012	4	599	6
156	4,112	2	982	3
625	18,658	2	2,635	3
2,500	101,526	4	10,389	6
10,000	457,225	5	50,694	11

Standard curve data is from a representative experiment.

Kit Size	SI2400	SI6000
1 plate	K15123B-1	K11123B-1
5 plates	K15123B-2	K11123B-2
20 plates	K15123B-3	K11123B-3
20 plates (Base)	K15123A-3	K11123A-3

Detection Limit (pg/mL)	rEPO	rVEGF
Average LLOD	2	31

LLOD is defined as 2.5x stdev above the background.