

# Measuring Immune Response and Cytokine Autoantibodies in Human Sepsis Samples

Robert Keith, Brooke Gilliam, Hong Luo, Harold Steiner, Yiwen Jan, Danielle Pepin, Qiang Xiao, and Wei Zheng  
MilliporeSigma, St. Louis, Missouri 63103



## Introduction

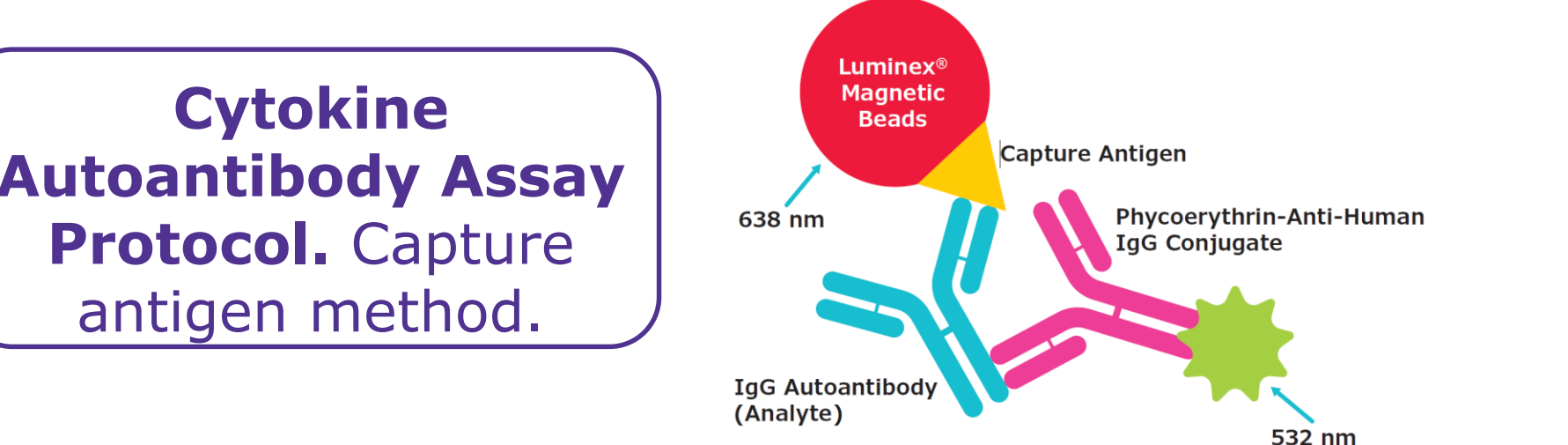
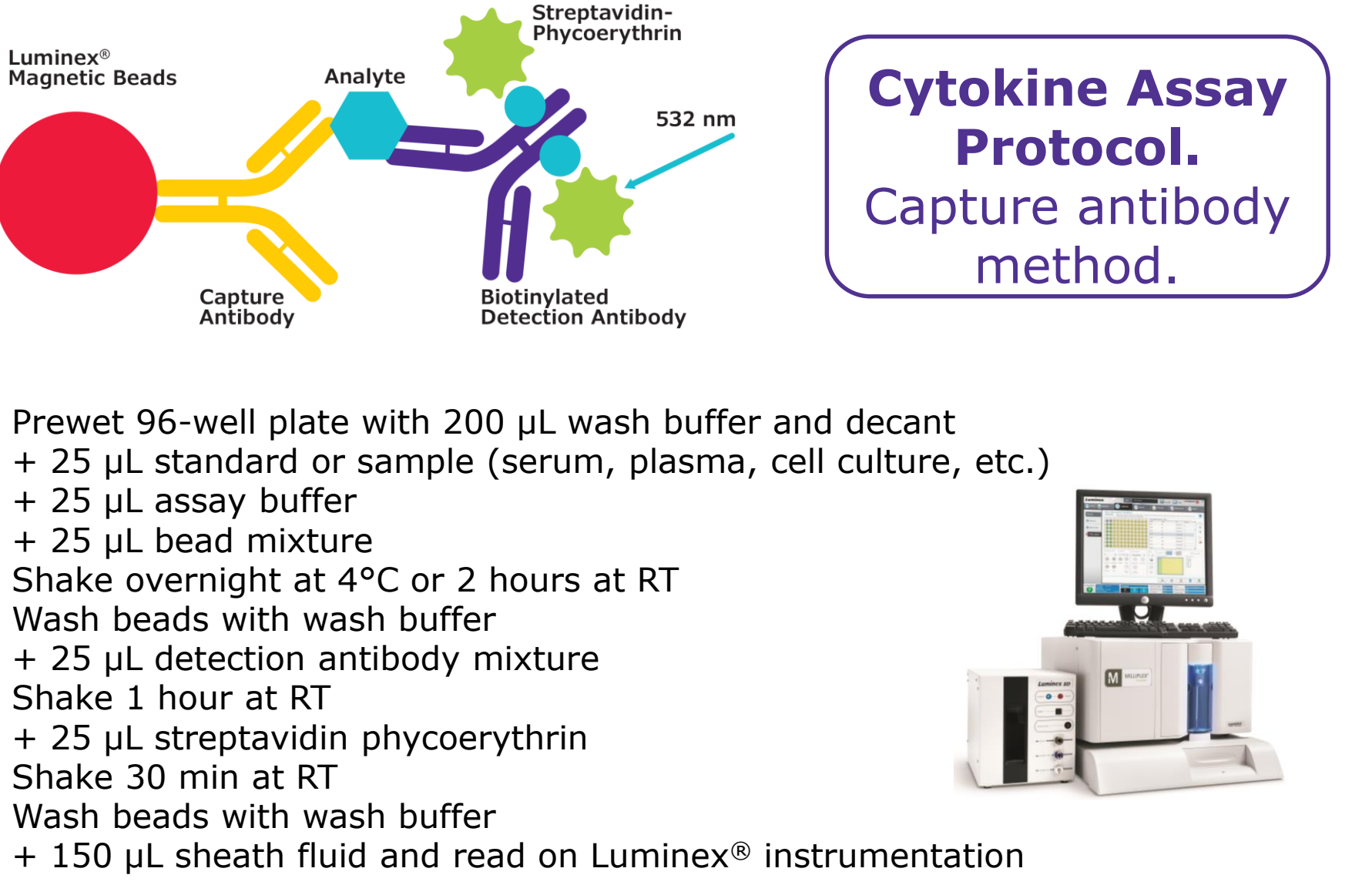
Cytokine, chemokine and growth factor research plays a significant role in achieving a deeper understanding of the immune system as well as many disease states such as inflammatory disease, allergic reactions, IBD, sepsis, and cancer. The ability to test for multiple factors simultaneously in a single sample is a valuable tool to researchers. MILLIPLEX® Human Cytokine/Chemokine/Growth Factor Panel A (Cat. No. [HCYTA-60K](#)) combines tests for 48 individual immune factors that previously have not been together in a single panel.

Anti-cytokine antibodies occur frequently and are present in healthy individuals and patients with acquired immunodeficiency and autoimmune diseases. Cytokines offering protection against microbes can be targeted by cytokine autoantibodies, leading to life-threatening infections. Measuring cytokine autoantibodies may be useful for disease monitoring and efficacy of treatment. We are developing a MILLIPLEX® Human Cytokine Autoantibody Panel (Cat. No. [HCYTAAB-17K](#)) consisting of 15 autoantibody immunoassays for use in serum and plasma samples. We measured expression of autoantibodies in serum and plasma from patients with SLE, RA, and sepsis along with healthy controls.

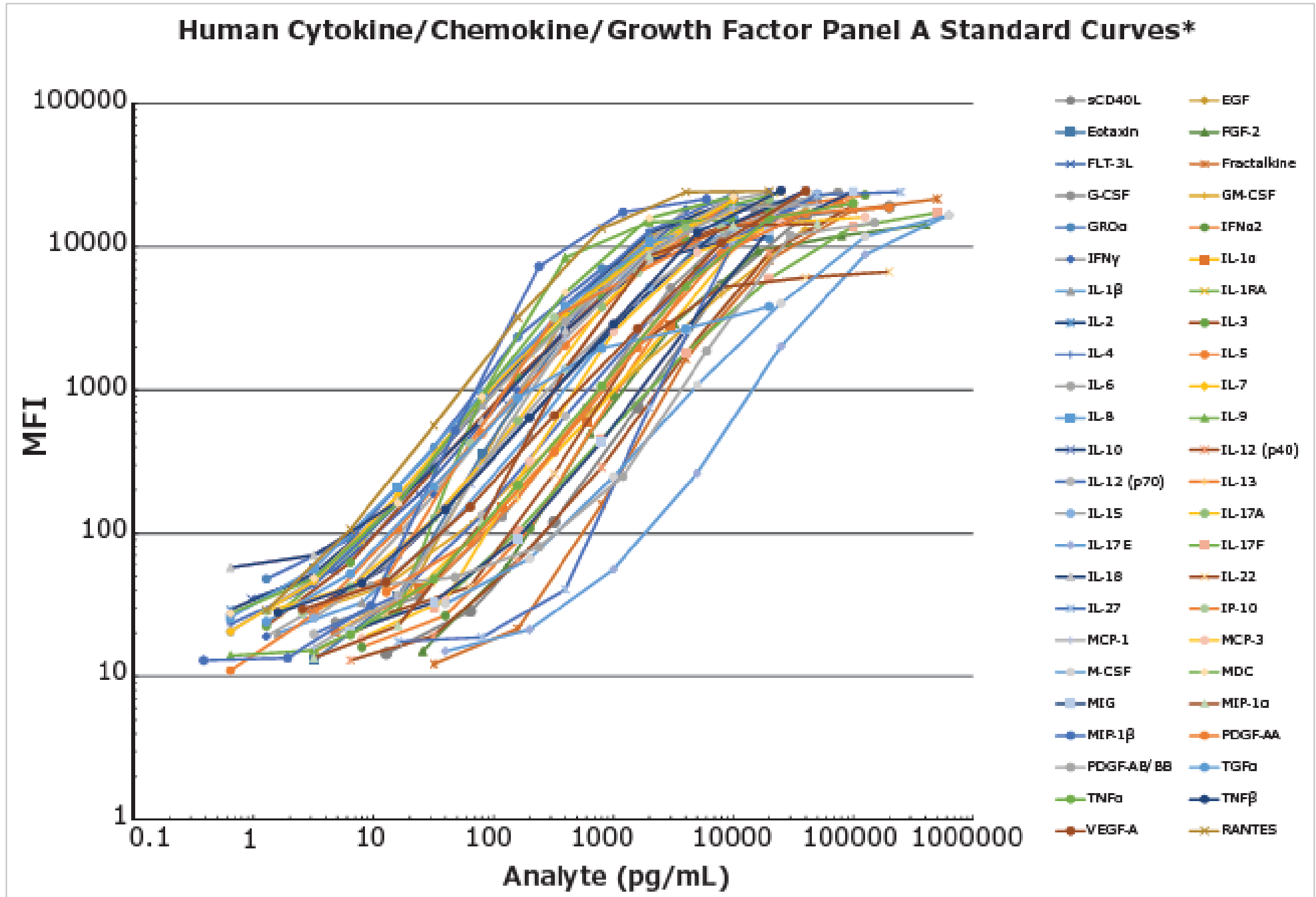
The use of these two panels allows researchers to gain information they may not have otherwise been able to access elucidating the relationship between cytokine levels and the corresponding autoantibody levels.

## Methods

**Microspheres.** We used magnetic microsphere beads from Luminex® Corp. Each set of beads is distinguished by different ratios of two internal dyes yielding a unique fluorescent signature to each bead set. Capture antibodies or antigens were coupled to the magnetic beads.



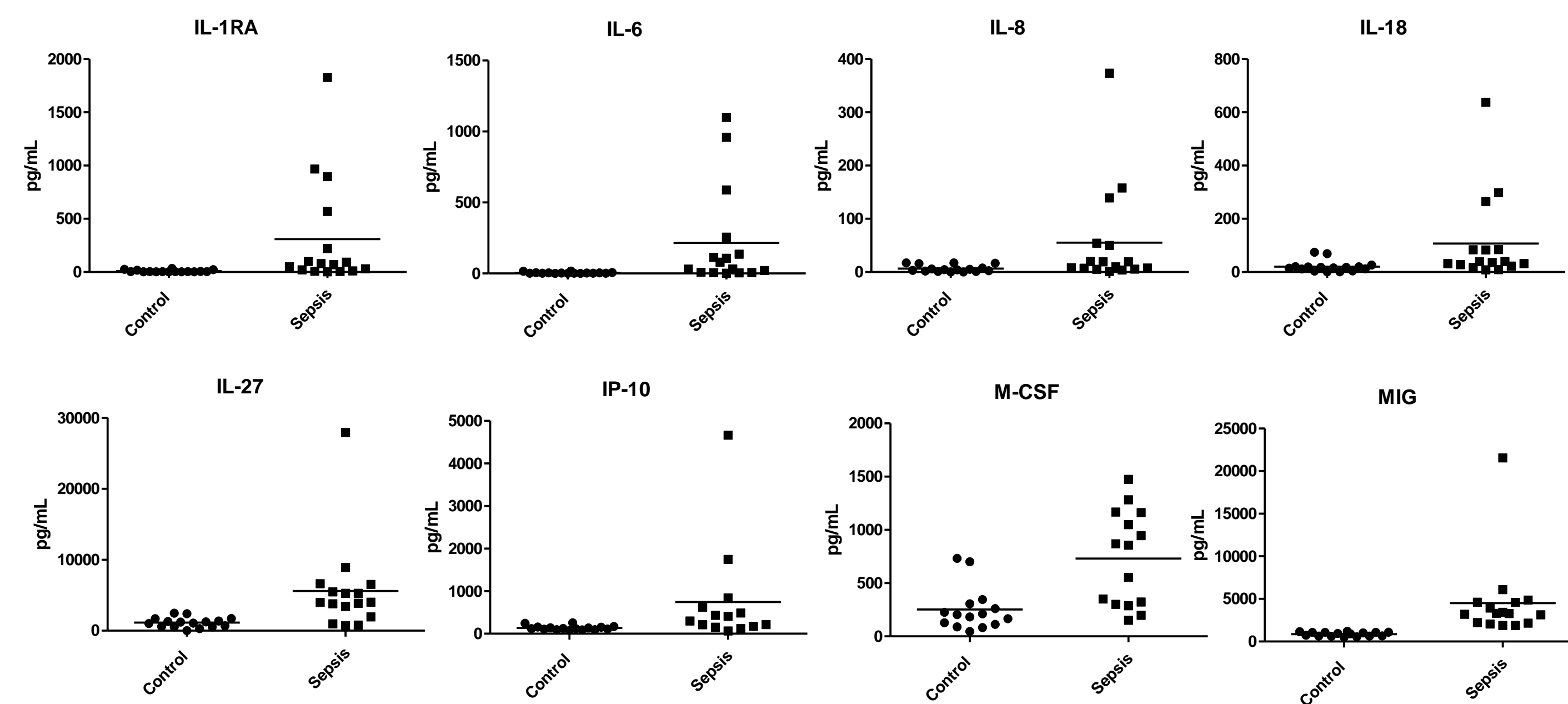
## 48-Plex Characteristics



\*Standard curves: Performed in MXHSM-A serum matrix except \*RANTES, which was in L-AB assay buffer. The kit was run using the [HCYTA-60K](#) overnight assay protocol.  
Sensitivity: Standard curves (n=9 separate tests) were run in serum matrix, except RANTES was tested in assay buffer.  
Precision: Two different standard controls were tested in serum matrix for intra-assay (n=8) and inter-assay (n=9) studies.  
Accuracy: Three different standard controls were spiked into serum matrix (n=6).

Analyte	Standard Curve Range (pg/mL)	Sensitivity (pg/mL)	Precision (%CV)		Accuracy (% Recovery) in Matrix
			Intra-assay	Inter-assay	
sCD40L	12.8-200,000	6.40	4.20	18.39	95
EGF	3.2-50,000	3.19	1.97	4.57	101
Eotaxin	3.2-50,000	3.07	3.93	5.27	96
FGF-2	29.6-400,000	22.40	4.73	4.33	106
Flt3 Ligand	0.96-15,000	0.85	3.11	4.66	95
Fractalkine	32-500,000	29.61	3.17	5.36	98
G-CSF	4.8-75,000	3.53	9.29	7.71	97
GM-CSF	2.98-40,000	1.60	5.82	4.78	100
GROα	1.28-20,000	1.06	3.96	3.93	99
IFNα2	8-125,000	6.43	4.35	4.39	98
IFNγ	1.28-20,000	0.86	10.10	12.74	94
IL-1α	4.8-75,000	2.22	5.00	5.11	98
IL-1β	1.6-25,000	0.52	3.85	5.92	92
IL-1RA	1.6-25,000	1.26	4.00	6.29	97
IL-2	0.64-10,000	0.27	4.02	4.82	95
IL-3	1.28-20,000	0.29	3.26	5.62	100
IL-4	0.64-10,000	0.20	4.55	5.94	95
IL-5	0.64-10,000	0.17	3.99	5.42	96
IL-6	0.64-10,000	0.14	5.06	6.13	96
IL-7	0.64-10,000	0.14	4.85	4.61	97
IL-8	0.64-10,000	0.52	3.07	5.05	96
IL-9	0.64-10,000	2.82	3.47	5.13	101
IL-10	2.56-40,000	0.88	3.91	5.88	98
IL-12 (p40)	6.4-100,000	3.19	3.35	4.68	95
IL-12 (p70)	3.2-50,000	0.88	3.53	5.99	96
IL-13	6.4-100,000	2.42	11.67	7.40	95
IL-15	3.2-50,000	0.71	4.91	5.69	96
IL-17A	1.28-20,000	0.70	4.72	5.14	93
IL-17E/IL-25	40-625,000	21.25	5.03	5.45	93
IL-17F	32-500,000	28.56	4.63	6.08	95
IL-18	0.64-10,000	0.53	4.38	4.68	97
IL-22	12.8-200,000	12.62	4.54	6.61	96
IL-27	16-250,000	47.50	7.24	4.82	96
IP-10	2.56-40,000	2.13	4.14	9.11	99
MCP-1	3.2-50,000	3.07	3.91	4.25	101
MCP-3	8-125,000	8.58	3.51	4.76	104
GM-CSF	40-625,000	31.94	4.38	5.04	97
MDC	0.64-10,000	0.42	3.47	4.27	96
MIG	6.4-100,000	3.97	6.75	11.07	95
MIP-1α	3.2-50,000	3.72	2.92	3.12	104
MIP-1β	0.38-6,000	0.34	3.56	4.73	99
PDGF-AA	12.8-200,000	10.31	5.88	5.98	102
PDGF-AB/BB	9.6-150,000	16.05	4.71	5.92	91
RANTES*	1.28-20,000	1.59	6.36	7.29	93
TGFα	1.28-20,000	0.98	3.80	5.39	102
TNFα	6.4-100,000	5.30	4.36	5.77	101
TNFβ	1.6-25,000	0.77	3.50	7.12	99
VEGF-A	2.56-40,000	0.96	4.94	5.73	99

## 48-Plex Sepsis vs. Normal Results



Healthy control serum/plasma samples (obtained from BioIVT) and sepsis patient serum/plasma samples (obtained from BioIVT, Discovery, and BioChemed) were tested neat (25 µL/well) in the [HCYTA-60K](#) panel.  
Healthy control serum/plasma samples, N=20. Sepsis serum/plasma samples, N=16.

Serum/Plasma Sample Testing: Serum/Plasma samples were tested neat (25 µL/well) except for RANTES for which samples were diluted 1:100 in assay buffer.

Normal Serum/Plasma Sample Summary Data (pg/mL)														
N	sCD40L	EGF	Eotaxin	FGF-2	FLT-3L	Fractalkine	G-CSF	GM-CSF	GROα	IFNα2	IFNγ	IL-1α	IL-1β	IL-1RA
average	2273.5	93.8	85.1	151.8	24.3	436.8	120.9	31.8	31.4	77.4	34.5	39.8	20.0	20.0
min	29.4	6.1	4.2	24.4	3.4	76.0	5.7	1.2	4.5	10.1	1.2	1.9	20.0	20.0
max	9505.0	353.7	126.8	499.2	81.6	1583.0	646.5	144.1	60.1	250.6	120.8	189.3	20.0	20.0
% detectable	100.0	85.0	100.0	70.0	100.0	75.0	90.0	70.0	80.0	65.0	85.0	75.0	85.0	75.0
Sepsis Serum/Plasma Sample Summary Data (pg/mL)														
N	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0
average	1886.9	63.2	70.3	68.6	28.9	214.8	1250.4	13.5	62.2	24.4	29.4	22.6	20.0	20.0
min	24.0	17.0	22.0	12.5	91.0	17.0	23.0	22.5	46.0	14.0	18.0	17.0	20.0	20.0
max	7097.4	196.5	137.2	138.0	50.9	446.5	11850.0	55.8	222.5	88.5	79.1	78.5	20.0	20.0
Normal Serum/Plasma Sample Summary Data (pg/mL)														
N	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0
average	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0	20.0
min	0.7	1.7	0.4	0.6	0.2	0.3	0.5	0.4	1.0	2.5	0.9	6.5	20.0	20.0
max	102.6	34.6	30.3	3.3	20.8	33.3	17.1	12.9	17.2	55.5	167.2	744.6	20.0	20.0
% detectable	85.0	100.0	85.0	45.0	90.0	100.0	95.0	100.0	90.0	100.0	95.0	100.0	85.0	100.0
Sepsis Serum/Plasma Sample Summary Data (pg/mL)														
N	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0
average	9.5	309.3	1.8	2.0	1.9	4.6	215.4	9.1	55.1	14.8	162.8	96.3	16.0	16.0
min	1.0	3.1	0.4	0.4	0.3	0.8	1.5	2.0	1.2	4.9	1.2	35.8	16.0	16.0
max	41.4	1828.2	10.7	5.9	7.5	28.6	1099.0	28.3	373.5	89.6	1796.5	247.1	16.0	16.0

Normal Serum/Plasma Sample Summary Data (pg/mL)														
N	IL-12 (p70)	IL-13	IL-15	IL-17A*	IL-17E	IL-17F	IL-18	IL-22	IL-27	IP-10	MCP-1	MCP-3		
average	33.9	72.2	17.1	27.6	1205.7	2630.4	20.4	331.2	1204.3	136.3	256.7	103.9		
min	2.1	3.0	3.0	0.8	52.5	1718.0	0.8	40.8	287.5	76.2	42.6	11.1		
max	119.0	451.3	101.0	56.2	6145.6	4313.0	74.0	863.2	2472.3	254.5	756.5	405.1		
% detectable	65.0	13.0	100.0	25.0	100.0	20.0	100.0	55.0	95.0	100.0	100.0	85.0		
Sepsis Serum/Plasma Sample Summary Data (pg/mL)														
N	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0		
average	6.1	23.9	17.3	5.0	487.0	212.7	107.1	75.0	5587.5	1663.7	573.2	31.3		
min	1.6	8.3	5.4	0.8	25.0	46.5	8.0	15.6	734.0	65.5	107.4	11.1		
max	31.2	70.2	40.3	19.4	2921.9	3079.0	637.6	218.2	27900.0	14532.0	2041.3	64.8		
Normal Serum/Plasma Sample Summary Data (pg/mL)														
N	M-CSF	MDC	MIG	MIP-1α	MIP-1β	PDGF-AA	PDGF-AB/RANTES	TGFα	Tnfα	TNFβ	VEGF-A			
average	20.0	20.0	20.0	20.0	20.0	20.0	20.0	8.0	20.0	20.0	20.0	20.0		
min	25.8	842.9	869.7	393.2	22.9	2272.4	15919.9	6983.5	118.5	82.4	72.0	130.6		
max	732.5	1592.0	1205.6	1434.5	39.7	7147.8	32160.0	129934.0	755.2	425.0	352.8	379.9		
% detectable	95.0	100.0	100.0	85.0	100.0	100.0	100.0	80.0	100.0	100.0	75.0	100.0		
Sepsis Serum/Plasma Sample Summary Data (pg/mL)														
N	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0	16.0		
average	731.3	729.0	4512.0	24.9	47.8	3285.0	20776.9	116456.5	24.0	61.1	8.1	173.1		
min	150.6	328.5	1881.5	5.4	33.5	1480.0	10630.4	26792.0	2.2	10.6	1.2	37.3		
max	1474.4	2802.0	21550.0	87.9	155.2	5573.2	47940.0	293916.0	137.6	293.5	32.8	489.9		

## Cytokine Autoantibody Results

Fifteen sepsis serum/plasma samples were