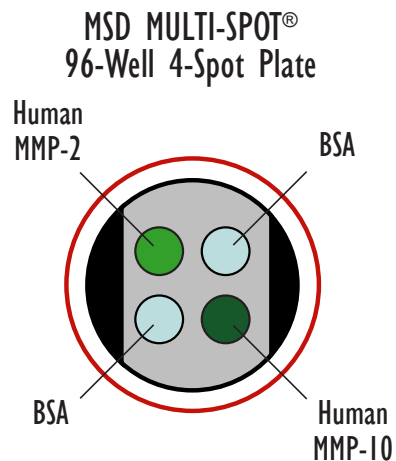
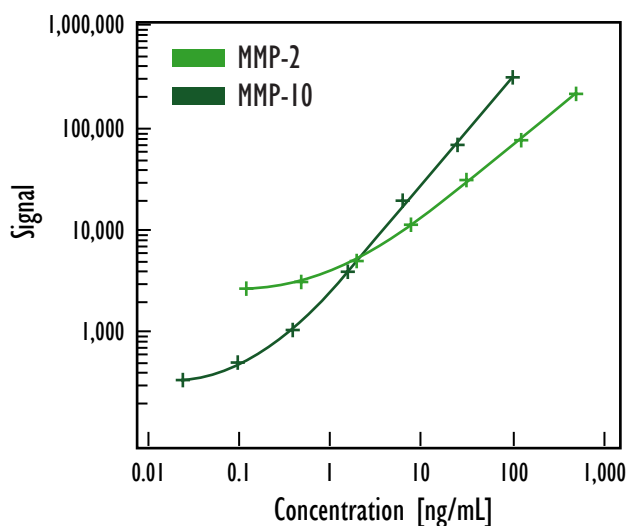


# MULTI-SPOT<sup>®</sup> Human Matrix Metalloprotease Ultrasensitive Assay

## Detection of MMP-2 and MMP-10 in Human Serum Samples



### Endogenous Levels (ng/mL)

Matrix	MMP-2	MMP-10
Serum	46.0	0.7
Heparin Plasma	36.9	1.1

Endogenous levels are based on measurements of pooled Human Serum and Heparin Plasma samples.

### Dilutional Linearity

Dilutional Linearity	Dilution Factor	% Recovery MMP-2	% Recovery MMP-10
Serum	1/4	118	98
	1/8	110	94
Heparin Plasma	1/4	136	92
	1/8	147	102

### Detection Limits

Detection Limit (ng/mL)	MMP-2	MMP-10
LLOD	1.4	1.4

Signal (Serum Protocol)	
Concentration (ng/mL)	MMP-2
500	233,968
125	71,075
31.25	31,578
7.8	11,153
2.0	4,655
0.5	2,903
0.12	2,698
0	2,190

Signal (Serum Protocol)	
Concentration (ng/mL)	MMP-10
100	276,042
25	63,813
6.25	18,397
1.6	3,791
0.4	1,128
0.1	495
0.02	347
0	269

### Recoveries

% Spike Recovery	Spike (ng/mL)	MMP-2	MMP-10
Serum	125	82	105
	31.3	104	144
	7.8	110	101
Heparin Plasma	125	77	105
	31.3	89	88
	7.8	96	100

Serum and plasma samples spiked with MMP-2 and MMP-10 were run undiluted.

Kit Size	SI2400	SI6000
1 plate	KI5033C-1	KI1033C-1
5 plates	KI5033C-2	KI1033C-2
20 plates	KI5033C-4	KI1033C-4
20 plates (Base)	KI5033A-1	KI1033A-1

# Multiplexing MMPs

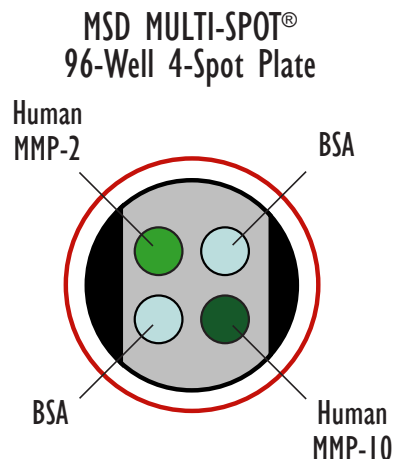
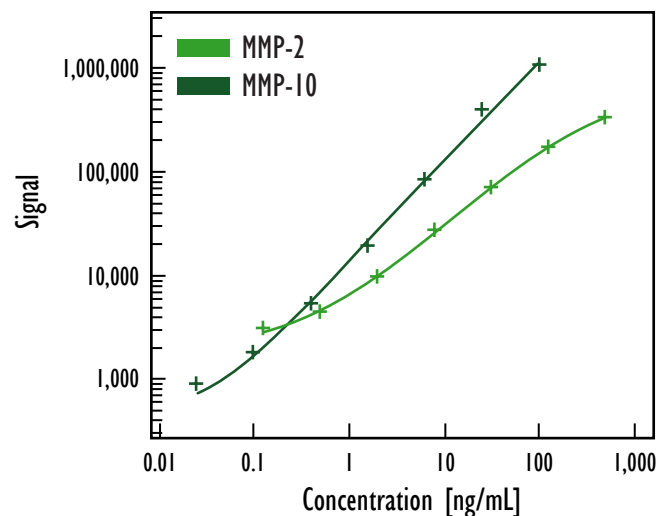
## Cross-Reactivity of MSD MMP Assays Using the Serum Protocol and the Human MMP 2-plex (2,10) Ultrasensitive Kit

MMP calibrators were assayed individually using the Human MMP 2-plex Ultrasensitive Kit.

Calibrator	Concentration (ng/mL)	Counts on Spot		% Cross-reactivity	
		MMP-2	MMP-10	MMP-2	MMP-10
MMP-2	0	2,375	260	100.00%	0.01%
	100	40,946	324		
MMP-10	0	2,375	260	2.96%	100.00%
	100	3,516	503,964		
MMP-1	0	2,375	260	-0.33%	0.00%
	100	2,246	262		
MMP-3	0	2,375	260	-0.18%	0.00%
	100	2,307	264		
MMP-9	0	2,375	260	-0.27%	0.00%
	100	2,269	281		
MMP-7	0	2,375	260	-0.02%	0.01%
	100	2,367	287		

# MULTI-SPOT<sup>®</sup> Human Matrix Metalloprotease Ultrasensitive Assay

## Detection of MMP-2 and MMP-10 in Tissue Culture Samples



### Detection Limits

Detection Limit (ng/mL)	MMP-2	MMP-10
LLOD	0.048	0.003

Signal (TCM Protocol)	
Concentration (ng/mL)	MMP-2
500	337,100
125	171,631
31.25	80,097
7.8	29,518
2.0	8,638
0.5	4,261
0.12	3,151
0	1,815

Signal (TCM Protocol)	
Concentration (ng/mL)	MMP-10
100	1,093,798
25	409,604
6.25	90,368
1.6	19,741
0.4	5,583
0.1	1,771
0.02	962
0	343

Kit Size	SI2400	SI6000
1 plate	KI5033C-1	KI1033C-1
5 plates	KI5033C-2	KI1033C-2
20 plates	KI5033C-4	KI1033C-4
20 plates (Base)	KI5033A-1	KI1033A-1