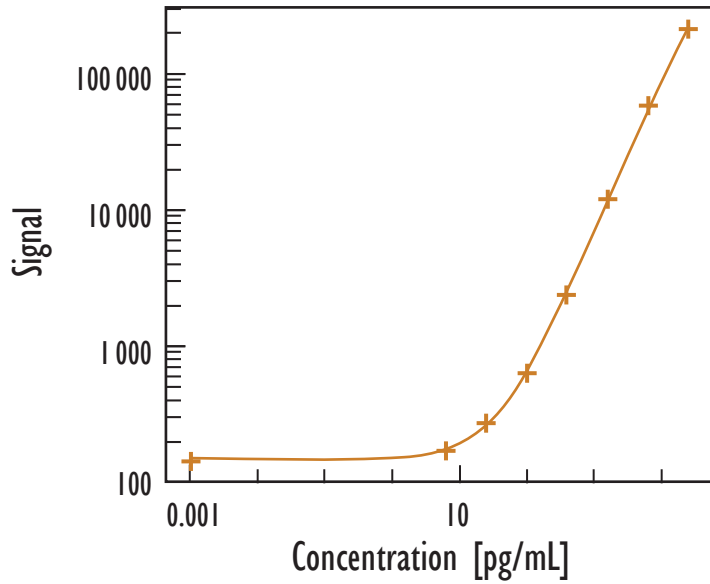


# MULTI-ARRAY<sup>®</sup> Human Cardiac Troponin T Assay

## Detection of Cardiac Troponin T in Human Serum and Plasma Samples



cTnT		
Concentration (pg/mL)	Mean Signal	%CV
0	144	10.2
2.4	174	6.0
9.8	277	9.1
39	643	6.4
156	2425	5.5
625	11981	3.0
2500	58653	4.8
10000	211107	2.7

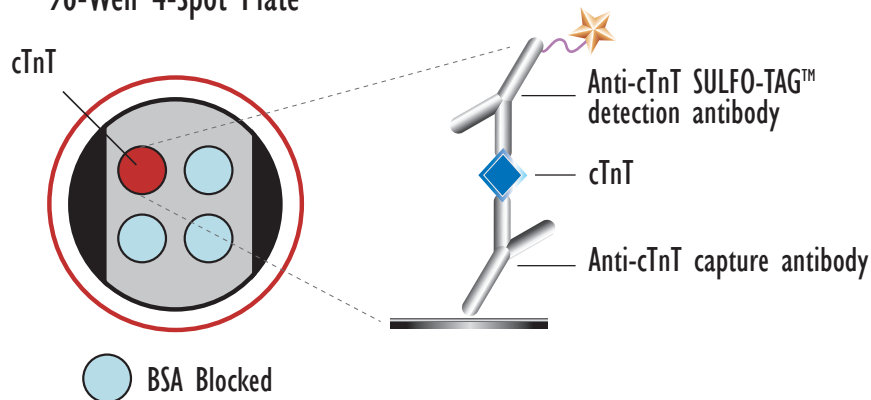
Standard curve data is from a representative experiment

Avg % CV is the average of CV's from calibrator levels above LLOD

cTnT LLOD (pg/mL)	7.2
-------------------	-----

LLOD (Lower Limit of Detection) is defined as 2.5x stdev above the background

MSD MULTI-SPOT<sup>®</sup>  
96-Well 4-Spot Plate



Kit Size	Catalog Number
1 plate	K151EFC-1
5 plates	K151EFC-2
20 plates	K151EFC-3
20 plates (Base)	K151EFA-3

# MULTI-ARRAY<sup>®</sup> Human Cardiac Troponin T Assay

## Detection of Cardiac Troponin T in Human Serum and Plasma Samples

### Recovery of Dilution Linearity

	Dilution Factor	Troponin T
Serum	1:2	96
	1:4	98
	1:8	107

- Samples were spiked with 4ng/mL Troponin T and diluted in calibrator diluent followed by subsequent dilution
- $\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$
- Values presented are averages across three pooled samples

### Endogenous Analyte Levels in Samples

	Troponin T (ng/mL)
N	20
Mean	< 0.005
Median	< 0.005
Range	< 0.005 - 0.006

- 20 normal human serum samples
- Average CVs for measured samples was less than 7%

### Recovery of Spiked Calibrator

	Spike Level (ng/mL)	Troponin T
Serum	0.5	113
	1.0	103
	2.0	109
EDTA Plasma	0.5	118
	1.0	107
	2.0	114
Heparin Plasma	0.5	114
	1.0	97
	2.0	109

- Measured analyte spiked into human samples
- $\% \text{ recovery} = \frac{(\text{measured value} * 100)}{\text{expected value}}$