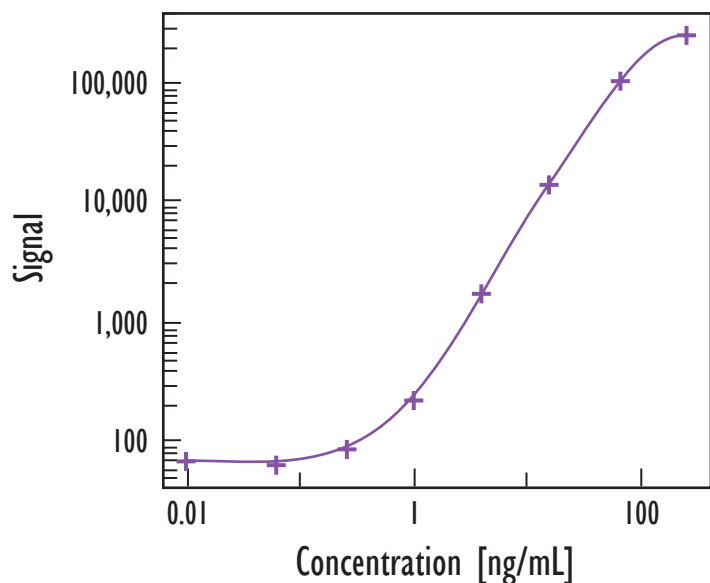


MULTI-ARRAY[®] Rat Cardiac Troponin ITC Complex Assay

Detection of Cardiac Troponin ITC Complex



ITC	
Concentration (ng/mL)	Mean Signal
0	64
0.06	60
0.24	83
0.98	204
3.91	1,652
15.6	13,750
62.5	103,339
250	256,861

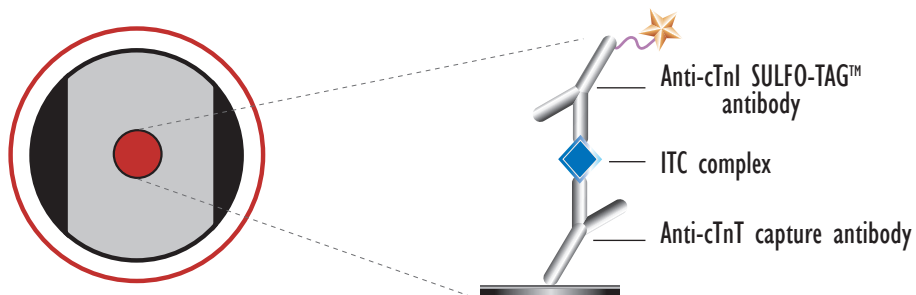
ITC LLOD (ng/mL)	0.42
------------------	------

Standard curve data is from a representative experiment

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

- The cardiac ITC complex assay does not cross-react with cardiac Tnl or cardiac TnT monomers.
- The capture antibody is cardiac TnT-specific, and the detection antibody is cardiac Tnl specific.

MSD MULTI-SPOT[®]
96-Well Small Spot Plate



Kit Size	Catalog Number
1 plate	K153HRC-1
5 plates	K153HRC-2
20 plates	K153HRC-3
20 plates (Base)	K153HRA-3

MULTI-ARRAY[®] Rat Cardiac Troponin ITC Complex Assay

Detection of Cardiac Troponin ITC Complex

Spike Recovery

- Normal, pooled rat serum was spiked with calibrators and assayed after 2-fold dilution into assay diluent

$$\% \text{ recovery} = \frac{(\text{measured value} * 100)}{\text{expected value}}$$

	Spike Level (ng/mL)	cITC
Serum	1.25	138%
	5	126%

Dilution Linearity

- Predicted value in calculation is based on measurement at 2-fold dilution
- Serum tested was pooled from treated rats

$$\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$$

	Dilution Factor	cITC
Serum	1/4	87%
	1/8	87%

Endogenous Levels

- Serum samples from normal and treated rats were tested at 2-fold dilution

	ng/mL	cITC
Serum	Normal	< LLOD
	Treated	3.6