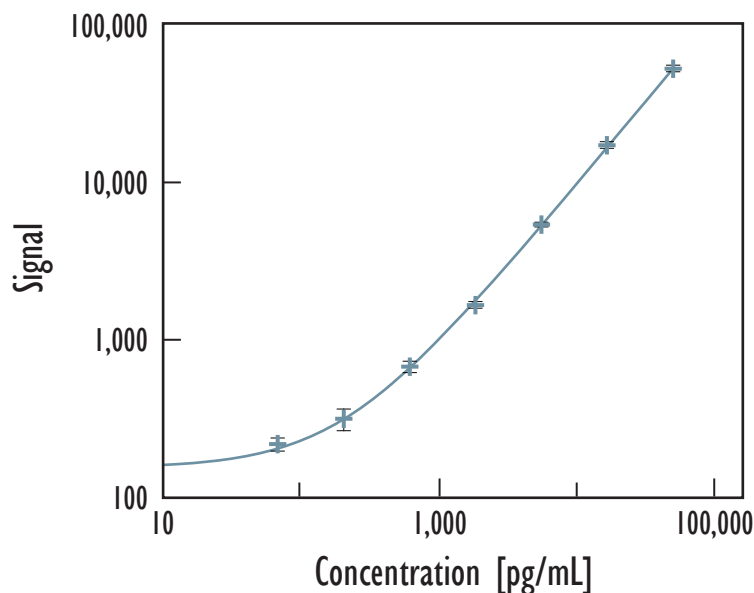


MULTI-ARRAY[®] Human Myeloperoxidase (MPO)

Detection of Myeloperoxidase in Human Serum and Plasma Samples



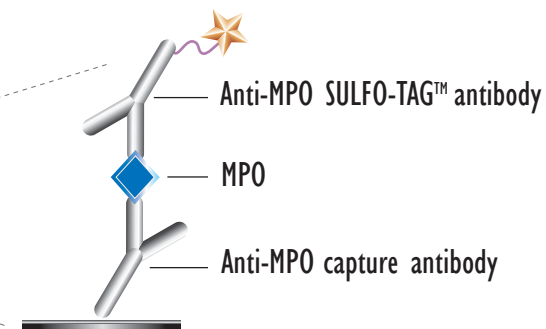
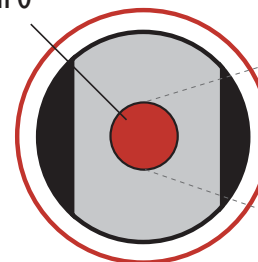
Concentration (pg/mL)	Signal
0	153
69	207
206	289
617	559
1,852	1,400
5,556	4,164
16,667	13,426
50,000	46,380

Above is representative calibration curve data

1:20 dilution of serum and plasma samples is recommended

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

MPO



Detection Limit (pg/mL)	MPO
LLOD	110

LLOD is defined as 2.5x stdev above the background

Kit Size	Catalog Number
1 plate	K151EEC-1
5 plates	K151EEC-2
20 plates	K151EEC-3

MULTI-ARRAY[®] Human Myeloperoxidase (MPO)

Detection of Myeloperoxidase in Human Serum and Plasma Samples

Spike Recovery

- Four serum samples were spiked with known amounts of MPO at defined levels and measured. The average spike recovery is tabulated.

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

% Recovery of Spiked Calibrator				
Spike Level (ng/mL)	Serum Samples			
	1	2	3	4
50	99	96	96	115
250	99	94	99	96
500	96	105	106	104

Dilutional Linearity

- Serum samples from three individuals each were diluted with the appropriate calibrator diluents. Recoveries were measured at three dilution levels – the average recovery at each dilution level is tabulated.

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

% Recovery of Dilution Linearity			
Dilution Factor	Serum Samples		
	1	2	3
1/2	98	100	90
1/4	102	103	97
1/8	106	94	89

Endogenous Levels in Human Samples

- Twenty serum and plasma (Sodium EDTA or Lithium Heparin) samples from healthy donors, approximately half female and half male were measured to obtain typical endogenous levels.
- Lithium Heparin and Sodium EDTA were found not to interfere with the MPO assay up to 50 U/mL and 50 mg/mL, respectively.

ng/mL	Serum	EDTA Plasma	Heparin Plasma
Mean	64.2	41.4	103.7
Median	57.4	32.3	98.5
Range	19.5 - 149.6	12.7 - 126.6	24.5 - 196.6